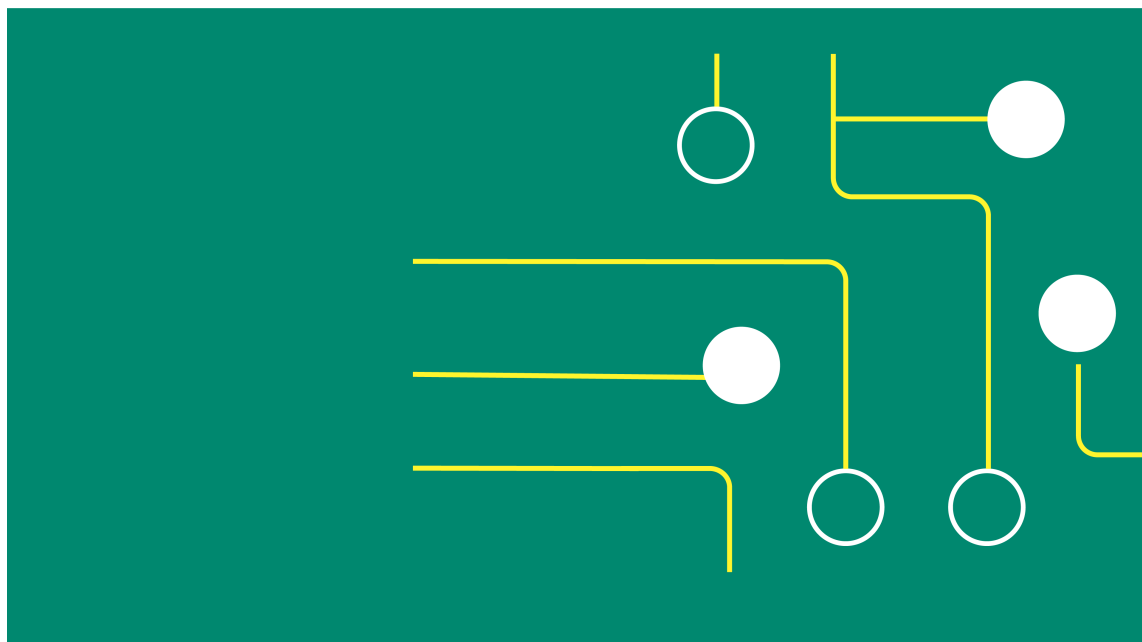




# UNICORN™ software

## Compatibility Matrix



# Table of Contents

<b>1</b>	<b>UNICORN 7.x compatibility .....</b>	<b>3</b>
1.1	Supported operating systems .....	4
1.2	Supported SQL Server and Databases .....	5
1.3	OPC Server compatibility .....	8
1.4	Instrument configurations compatibility .....	9
1.4.1	<i>ÄKTA avant 150</i> .....	10
1.4.2	<i>ÄKTA avant 25</i> .....	11
1.4.3	<i>ÄKTA pure 150</i> .....	12
1.4.4	<i>ÄKTA pure 25</i> .....	13
1.4.5	<i>Several ÄKTA systems, ReadyToProcess, and WAVE</i> .....	14
1.4.6	<i>Legacy ÄKTA systems</i> .....	16
<b>2</b>	<b>UNICORN 6.x compatibility .....</b>	<b>17</b>
2.1	Supported operating systems .....	18
2.2	Supported SQL Server and databases .....	19
2.3	Instrument configurations compatibility .....	20
2.3.1	<i>ÄKTA avant 150</i> .....	21
2.3.2	<i>ÄKTA avant 25</i> .....	22
2.3.3	<i>ÄKTA pure 150</i> .....	23
2.3.4	<i>ÄKTA pure 25</i> .....	24
2.3.5	<i>Several ÄKTA systems, ReadyToProcess WAVE, and UniFlux</i> .....	25
2.3.6	<i>Legacy ÄKTA systems</i> .....	26
<b>3</b>	<b>Support for RTU and RTU Win10 Image .....</b>	<b>27</b>
<b>4</b>	<b>Modifications history .....</b>	<b>28</b>

# 1 UNICORN 7.x compatibility

## In this chapter

<b>Section</b>		<b>See page</b>
1.1	Supported operating systems	4
1.2	Supported SQL Server and Databases	5
1.3	OPC Server compatibility	8
1.4	Instrument configurations compatibility	9

## 1.1 Supported operating systems

UNICORN™ version	Windows Operating System					
	Win 7 SP1	Win 10 Pro 64-bit verified versions <sup>1</sup>	Win 10 Ent 64-bit verified versions <sup>1</sup>	Win 11 Pro 64-bit verified versions <sup>1</sup>	Win 11 Ent 64-bit verified versions <sup>1</sup>	OS Language
7.11 <sup>2,3</sup> , 7.11 SP1 <sup>2,3</sup>	N/A	22H2	22H2, LTSC 2019, LTSC 2021	22H2	22H2, 23H2	US/UK English
7.10 <sup>2</sup> , 7.10 AB <sup>2,3</sup> , 7.10 SP1 <sup>2,3</sup> , SP2 <sup>2,3</sup> , SP3 <sup>2,3</sup>	N/A	21H2	21H2, 2016LTSB, 2019LTSC	22H2	21H2	US/UK English
7.9 <sup>2</sup> , 7.9 SP2 <sup>2,3</sup>	N/A	21H2	21H2, 2016LTSB, 2019LTSC	22H2	21H2	US/UK English
7.8 <sup>2</sup> , 7.8 SP1 <sup>2,3</sup>	N/A	21H1	21H2, 2016LTSB, 2019LTSC	N/A	N/A	US/UK English
7.7 <sup>2</sup> , 7.7 SP1 <sup>2,3</sup>	N/A	20H2	21H2, 2016LTSB, 2019LTSC	N/A	N/A	US/UK English
7.6 SP1 <sup>2</sup> , SP2 <sup>2,3</sup>	N/A	20H2	20H2	N/A	N/A	US/UK English
7.6 <sup>2</sup>	N/A	1909	1909, 2016LTSB, 2019LTSC	N/A	N/A	US/UK English
7.5 SP1 <sup>2</sup>	N/A	20H2	1809, 1903	N/A	N/A	US/UK English
7.5 <sup>2</sup>	N/A	1809	1809, 1903	N/A	N/A	US/UK English
7.4 <sup>2</sup>	64-bit	1809	1709	N/A	N/A	US/UK English
7.3 SP2	64-bit	1809, 1909	2016 1607 LTSC	N/A	N/A	US English
7.3 SP1	64-bit	1803	2016 1607 LTSC	N/A	N/A	US English
7.3	64-bit	1709	2016 1607 LTSC	N/A	N/A	US English
7.2	32/64-bit	1511	2015, 2016 1607	N/A	N/A	US English
7.1 SP2	32/64-bit	20H2	-	N/A	N/A	US English
7.1 SP1	32/64-bit	1511	-	N/A	N/A	US English
7.1	32/64-bit	-	-	N/A	N/A	US English
7.0.2 SP3, SP4	32/64-bit	1809	-	N/A	N/A	US English
7.0.2 SP2	32/64-bit	1703	-	N/A	N/A	US English
7.0.2 SP1	32/64-bit	1511	-	N/A	N/A	US English
7.0.2	32/64-bit	-	-	N/A	N/A	US English
7.0.1, 7.0.1 SP1	32/64-bit	N/A	N/A	N/A	N/A	US English
7.0, 7.0 SP1	32/64-bit	N/A	N/A	N/A	N/A	US English

<sup>1</sup> UNICORN is verified on these versions but other versions may also work

<sup>2</sup> The .NET version used requires Win 10 1607 or later

<sup>3</sup> UNICORN versions released after June 2023 require the root certificate that is part of Windows since May 2020

## 1.2 Supported SQL Server and Databases

The default database version installed with UNICORN is written in **BOLD**.

UNICORN version	SQL Server Express Edition	Cloud-DB	Database version	SQL Server tested combination						
				2008 R2	2012 SP1	2014 SP1	2016 SP2	2017	2019	2022
7.11, 7.11 SP1	2014, 2017, 2019, 2022	Azure Managed Instance Azure SQL DB AWS RDS SQL Server	<b>1.8.6</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>
7.10 SP2, SP3	2014, 2017, 2019	Azure Managed Instance Azure SQL DB AWS RDS SQL Server	1.8.6, <b>1.8.5</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.10	2014, 2017, 2019	Azure Managed Instance Azure SQL DB AWS RDS SQL Server	<b>1.8.5</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.9 SP2	2014, 2017, 2019	Azure Managed Instance Azure SQL DB AWS RDS SQL Server	1.8.6, 1.8.5, <b>1.8.4</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.9	2014, 2017, 2019	Azure Managed Instance Azure SQL DB AWS RDS SQL Server	1.8.5, <b>1.8.4</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.8 SP1	2014, 2017, 2019	N/A	1.8.6, 1.8.5, 1.8.4, <b>1.8.3</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	

1 UNICORN 7.x compatibility  
 1.2 Supported SQL Server and Databases

UNICORN version	SQL Server Express Edition	Cloud-DB	Database version	SQL Server tested combination						
7.8	2014, 2017, 2019	N/A	1.8.5, 1.8.4, <b>1.8.3</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.7SP1	2014, 2017, 2019	N/A	1.8.6, 1.8.5, 1.8.4, 1.8.3, <b>1.8.2</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.7	2014, 2017, 2019	N/A	1.8.5, 1.8.4, 1.8.3, <b>1.8.2</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.6SP2	2014, 2017	N/A	1.8.6, 1.8.5, 1.8.4, 1.8.3, 1.8.2, <b>1.8.1</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.6, 7.6SP1	2014, 2017	N/A	1.8.5, 1.8.4, 1.8.3, 1.8.2, <b>1.8.1</b>				X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	
7.5SP1	2014, 2017	N/A	1.8.5, 1.8.4, 1.8.3, 1.8.2, 1.8.1, <b>1.8.0</b>		X	X	X	X		
7.5	2017	N/A	1.8.5, 1.8.4, 1.8.3, 1.8.2, 1.8.1, <b>1.8.0</b>		X	X	X	X		
7.4	2014	N/A	<b>1.7.5</b>		X	X	X	X		
7.3SP2	2014	N/A	1.8.1, 1.7.5, <b>1.7.4</b>		X	X	X	X		
7.3, 7.3SP1	2014	N/A	1.7.5, <b>1.7.4</b>		X	X	X	X		
7.2	2014	N/A	<b>1.7.3</b> to 1.7.5		X	X				
7.1, 7.1 SP1, SP2	2014	N/A	<b>1.7.2</b> to 1.7.5		X	X				
7.0.2SP2, SP3, SP4	2014	N/A	1.7.5, 1.7.4, <b>1.7.1</b>		X	X				
7.0.2SP1	2014	N/A	<b>1.7.1</b>		X	X				

UNICORN version	SQL Server Express Edition	Cloud-DB	Data-base version	SQL Server tested combination						
7.0.2	2014	N/A	<b>1.7.1</b>		X	X	X	X		
7.0.1, 7.0.1 SP1	2008 R2	N/A	1.7.1, <b>1.7.0</b>	X	X	X				
7.0 SP1	2008 R2	N/A	1.7.1, <b>1.7.0</b>	X	X	X				
7.0	2008 R2	N/A	1.7.1, <b>1.7.0</b>	X						

<sup>1</sup> The following links refer to Microsoft pages concerning dependencies between Operating Systems and SQL Servers:

- [Minimum SQL Server version requirements for Windows 10 and Windows Server 2016.](#)
- [SQL Server 2019: Hardware and software requirements.](#)
- [SQL Server 2016 and 2017: Hardware and software requirements.](#)

- Note:**
- *UNICORN Database SP1 replaces the UNICORN Database patch UNICORN 7.0.2-7.10 DB Patch 1 released earlier.*
  - *UNICORN Database SP1 is compatible with UNICORN Database versions 1.7.1 and later.*

## 1.3 OPC Server compatibility

OPC Server version	UNICORN version	OPC standards		
		A&E	DA	HDA
1.0	7.8	Yes	Yes	Yes
1.0 SP1	7.8 + 7.8 SP1 <sup>1</sup>	Yes	Yes	Yes
1.1	7.9	Yes	Yes	Yes
1.1 SP1	7.9 + 7.9 SP2 <sup>1</sup>	Yes	Yes	Yes
1.2	7.10	Yes	Yes	Yes
1.2.1	7.10 + 7.10 SP1 <sup>1</sup>	Yes	Yes	Yes
1.2.1 SP1	7.10 + 7.10 SP2 <sup>1</sup>	Yes	Yes	Yes
1.3	7.11	Yes	Yes	Yes

<sup>1</sup> Both need to be installed



## 1.4 Instrument configurations compatibility

### In this section

<b>Section</b>	<b>See page</b>
1.4.1 ÄKTA avant 150	10
1.4.2 ÄKTA avant 25	11
1.4.3 ÄKTA pure 150	12
1.4.4 ÄKTA pure 25	13
1.4.5 Several ÄKTA systems, ReadyToProcess, and WAVE	14
1.4.6 Legacy ÄKTA systems	16

## 1.4.1 ÄKTA avant 150

ÄKTA avant™ 150 IC <sup>1</sup>	UNICORN version													
	7.0	7.0.1	7.0.2	7.1	7.2	7.3	7.4	7.5	7.6	7.7	7.8	7.9	7.10	7.11
1.0.5.0	X	X	X	X										
1.1.2.0	X	X	X	X										
1.2.5.0	X	X	X	X										
1.3.1.0	X	X	X	X										
1.4.0.0	X	X	X	X										
1.5.1.0	X	X	X	X										
1.6.0.0	X	X	X	X										
2.0.2.1	X	X	X	X										
2.1.0.2	X	X	X	X										
2.2.0.0	X	X	X	X										
2.3.0.0	X	X	X	X										
2.4.0.4	X	X	X	X	X	X	X <sup>2</sup>	X	X	X	X			
2.5.0.0	X	X	X	X	X	X	X <sup>2</sup>	X	X	X	X			
2.6.0.3	X	X	X	X	X	X	X <sup>2</sup>	X	X	X	X			
2.7.0.5	X	X	X	X	X	X	X <sup>2</sup>	X	X	X	X			
3.0.0.8									X	X	X			
3.1.0.30									X	X	X			
3.2.0.6									X <sup>3</sup>	X <sup>3</sup>	X <sup>3</sup>	X <sup>3,4</sup>	X <sup>3</sup>	X

<sup>1</sup> Instrument Configuration

<sup>2</sup> Due to a defect in the implementation of the DoE functionality in UNICORN 7.4, it is recommended to operate the system with a later version of UNICORN.

<sup>3</sup> For IC systems using a Real-Time Unit (RTU), an update of the RTU Win 10 image is needed (for image versions 2.0 or earlier): RTU Win 10 image SP1 must be installed.

<sup>4</sup> It is not recommended to install UNICORN 7.9 due to a change in how method parameters are handled. If UNICORN 7.9 is already installed, it is not recommended to update ICs. For details, see the *UNICORN 7.9 Method Editor Information Letter*.

## 1.4.2 ÄKTA avant 25

ÄKTA avant 25 IC	UNICORN version													
	7.0	7.0.1	7.0.2	7.1	7.2	7.3	7.4	7.5	7.6	7.7	7.8	7.9	7.10	7.11
2.0.5.0	X	X	X	X										
2.1.2.0	X	X	X	X										
2.2.5.0	X	X	X	X										
2.3.1.0	X	X	X	X										
2.4.1.0	X	X	X	X										
2.5.0.0	X	X	X	X										
3.0.2.1	X	X	X	X										
3.1.0.3	X	X	X	X										
3.2.0.0	X	X	X	X										
3.3.0.0	X	X	X	X										
3.4.0.15	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
3.5.0.0	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
3.6.0.15	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
3.7.0.4	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
4.0.0.8									X	X	X			
4.1.0.30									X	X	X			
4.2.0.16									X <sup>2</sup>	X <sup>2</sup>	X <sup>2</sup>	X <sup>2,3</sup>	X <sup>2</sup>	X

<sup>1</sup> Due to a defect in the implementation of the DoE functionality in UNICORN 7.4, it is recommended to operate the system with a later version of UNICORN.

<sup>2</sup> For IC systems using a Real-Time Unit (RTU), an update of the RTU Win 10 image is needed (for image versions 2.0 or earlier): RTU Win 10 image SP1 must be installed.

<sup>3</sup> It is not recommended to install UNICORN 7.9 due to a change in how method parameters are handled. If UNICORN 7.9 is already installed, it is not recommended to update ICs. For details, see the *UNICORN 7.9 Method Editor Information Letter*.

## 1.4.3 ÄKTA pure 150

ÄKTA pure™ 150 IC	UNICORN version													
	7.0	7.0.1	7.0.2	7.1	7.2	7.3	7.4	7.5	7.6	7.7	7.8	7.9	7.10	7.11
1.0.0.1	X	X	X	X										
1.2.0.3	X	X	X	X										
1.3.0.1	X	X	X	X										
1.4.1.2	X	X	X	X										
1.5.0.0	X	X	X	X										
1.6.0.2	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
1.7.0.6	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
1.8.0.3	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
2.0.0.4									X	X	X			
2.1.0.13									X	X	X			
2.2.0.3									X	X	X			
2.3.0.7									X <sup>2</sup>	X <sup>2</sup>	X <sup>2</sup>	X <sup>2,3</sup>	X <sup>2</sup>	X

<sup>1</sup> Due to a defect in the implementation of the DoE functionality in UNICORN 7.4, it is recommended to operate the system with a later version of UNICORN.

<sup>2</sup> For IC systems using a Real-Time Unit (RTU), an update of the RTU Win 10 image is needed (for image versions 2.0 or earlier): RTU Win 10 image SP1 must be installed.

<sup>3</sup> It is not recommended to install UNICORN 7.9 due to a change in how method parameters are handled. If UNICORN 7.9 is already installed, it is not recommended to update ICs. For details, see the *UNICORN 7.9 Method Editor Information Letter*.

## 1.4.4 ÄKTA pure 25

ÄKTA pure 25 IC	UNICORN version													
	7.0	7.0.1	7.0.2	7.1	7.2	7.3	7.4	7.5	7.6	7.7	7.8	7.9	7.10	7.11
1.0.1.2	X	X	X	X										
1.1.4.0	X	X	X	X										
1.2.0.0	X	X	X	X										
1.4.0.3	X	X	X	X										
1.5.0.2	X	X	X	X										
1.6.1.2	X	X	X	X										
1.7.0.1	X	X	X	X										
1.8.0.6	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
1.9.0.11	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
1.10.0.6	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
1.12.0.6	X	X	X	X	X	X	X <sup>1</sup>	X	X	X	X			
1.13.0.63	X	X	X	X	X <sup>2</sup>	X <sup>2</sup>	X <sup>1,2</sup>	X <sup>2</sup>	X	X	X			
2.0.0.20									X	X	X			
2.1.0.1									X	X	X			
2.2.0.27									X	X	X			
2.3.0.3									X	X	X			
2.4.0.8									X <sup>3</sup>	X <sup>3</sup>	X <sup>3</sup>	X <sup>3,4</sup>	X <sup>3</sup>	X

<sup>1</sup> Due to a defect in the implementation of the DoE functionality in UNICORN 7.4, it is recommended to operate the system with a later version of UNICORN.

<sup>2</sup> This IC has a certificate incompatible with older RTUs. Therefore, the IC is only valid for systems without RTU.

<sup>3</sup> For IC systems using a Real-Time Unit (RTU), an update of the RTU Win 10 image is needed (for image versions 2.0 or earlier): RTU Win 10 image SP1 must be installed.

<sup>4</sup> It is not recommended to install UNICORN 7.9 due to a change in how method parameters are handled. If UNICORN 7.9 is already installed, it is not recommended to update ICs. For details, see the *UNICORN 7.9 Method Editor Information Letter*.

**Note:** ÄKTA pure micro is supported by Instrument Configuration ÄKTA pure 25 version 2.0.0.20 or later.

## 1.4.5 Several ÄKTA systems, ReadyToProcess, and WAVE

Instrument configuration	UNICORN version																
	7.0	7.0.1	7.0.1SP1	7.0.2	7.0.2SP1	7.1	7.1SP1	7.2	7.3	7.4	7.5	7.6	7.7	7.8	7.9	7.10	7.11
ÄKTA pilot™ 600 1.0.0.53									X	X	X	X	X	X			
ÄKTA pilot 600 1.1.0.6									X	X	X	X	X	X			
ÄKTA pilot 600 1.2.0.2												X	X	X			
ÄKTA pilot 600 1.4.0.1												X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	X <sup>1,2</sup>	X <sup>1</sup>	X
ÄKTA go™ 1.0.0.17										X	X	X	X	X			
ÄKTA go 2.0.0.44												X	X	X			
ÄKTA go 2.1.0.0												X	X	X			
ÄKTA go 2.2.0.1												X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>			
ÄKTA go 3.0.0.4												X <sup>1</sup>	X <sup>1</sup>	X <sup>1</sup>	X <sup>1,2</sup>	X <sup>1</sup>	X
ÄKTA process™ CFG																X	X
(ÄKTA process™) 29165862_AA_W230(1.0.0.0)	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>2</sup>	X	X
Custom bioprocess	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>2</sup>	X	X
(ÄKTA ready™) Ready_202_W210 1.0.2.0		X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>2</sup>	X	X
Ready_203_W220 1.0.0.0		X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>2</sup>	X	X
Ready_205_W230 1.0.0.0		X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>2</sup>	X	X
ÄKTA ready XL				X	X	X	X	X	X	X	X	X	X	X	X <sup>2</sup>	X	X
<b>(ReadyToProcess™) WAVE</b> 25/Xuri™ 1.0.3.0	X	X	X	X	X	X	X										
ReadyToProcess WAVE™ 25/ Xuri 2.0.4.0	X	X	X	X	X	X	X				X	X	X	X	X <sup>2</sup>	X	X
WAVE 25/Xuri 2.0.7.0	X	X	X	X	X	X	X	X <sup>3</sup>	X <sup>3</sup>		X	X	X	X	X <sup>2</sup>	X	X
WAVE 25/Xuri 2.0.8.0	X	X	X	X	X	X	X	X <sup>3</sup>	X <sup>3</sup>		X	X	X	X	X <sup>2</sup>	X	X
WAVE 25/Xuri 2.0.9.0	X	X	X	X	X	X	X	X <sup>3</sup>	X <sup>3</sup>		X	X	X	X	X <sup>2</sup>	X	X
WAVE 25/Xuri 3.0.0.1													X	X	X <sup>2</sup>	X	X
WAVE 25/Xuri 3.1.0.12													X	X	X <sup>2</sup>	X	X
(ÄKTA readyflux™) ready- flux_1_0_5						X		X	X	X	X	X	X	X	X <sup>2</sup>	X	X
(ÄKTA readyflux) ready- flux_1_0_6								X	X	X	X	X	X	X	X <sup>2</sup>	X	X
(ÄKTA readyflux) ready- flux_2_0_0_3												X	X	X	X <sup>2</sup>	X	X
ÄKTA readyflux XL												X	X	X	X <sup>2</sup>	X	X
ÄKTA pcc™ 2.0.1.5				X				X	X	X	X	X	X	X	X <sup>2</sup>	X	X
ÄKTA pcc 2.1.0.11				X				X	X	X	X	X	X	X	X <sup>2</sup>	X	X
ÄKTA pcc 2.2.0.12				X				X	X	X	X	X	X	X	X <sup>2</sup>	X	X

# 1 UNICORN 7.x compatibility

## 1.4 Instrument configurations compatibility

### 1.4.5 Several ÄKTA systems, ReadyToProcess, and WAVE

Instrument configuration	UNICORN version																		
ÄKTA pcc 2.3.0.2															X	X	X <sup>2</sup>	X	X
ÄKTA pcc 2.4.0.10															X <sup>1</sup>	X <sup>1</sup>	X <sup>1,2</sup>	X <sup>1</sup>	X
ÄKTA oligosynt™ 1.0.0.44																X	X <sup>2</sup>	X	X
ÄKTA oligosynt 1.1.0.4																X	X <sup>2</sup>	X	X
ÄKTA oligosynt 1.2.0.6																X	X <sup>2</sup>	X	X
ÄKTA oligosynt 1.3.0.0																X <sup>1</sup>	X <sup>1,2</sup>	X <sup>1</sup>	X
OligoProcess™																X	X <sup>2</sup>	X	X
OligoPilot™																X	X <sup>2</sup>	X	X

<sup>1</sup> For IC systems using a Real-Time Unit (RTU), an update of the RTU Win 10 image is needed (for image versions 2.0 or earlier): RTU Win 10 image SP1 must be installed.

<sup>2</sup> It is not recommended to install UNICORN 7.9 due to a change in how method parameters are handled. If UNICORN 7.9 is already installed, it is not recommended to update ICs. For details, see the *UNICORN 7.9 Method Editor Information Letter*.

<sup>3</sup> Because of incompatibility between Griffin and UNICORN 7.2 and 7.3, these combinations are not recommended.

## 1.4.6 Legacy ÄKTA systems

Instrument configuration	UNICORN version																
	7.0	7.0.1	7.0.1SP1	7.0.2	7.0.2SP1	7.1	7.1SP1	7.2	7.3	7.4	7.5	7.6	7.7	7.8	7.9	7.10	7.11
(ÄKTA crossflow™) cross-flow_111_W105 1.0.0.17								X	X	X	X	X	X	X	X <sup>1</sup>	X	
UniFlux™	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	X
(ÄKTApilot) PiloX_133_W253_1.0.0.2	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
<b>(ÄKTAexplorer)</b> E10_414_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
E10F414_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
E100_444_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
E100F444_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
<b>(ÄKTApurifier/Basic)</b> P10F414_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
P10_415_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
P100_415_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
P100F415_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
<b>(ÄKTApurifier UPC)</b> P10_UPC_414_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
P10F_UPC_414_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
P100_UPC_414_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	
P100F_UPC_414_1.0.0.0	X	X	X	X	X	X	X	X	X	X	X	X	X	X	X <sup>1</sup>	X	

<sup>1</sup> It is not recommended to install UNICORN 7.9 due to a change in how method parameters are handled. If UNICORN 7.9 is already installed, it is not recommended to update ICs. For details, see the *UNICORN 7.9 Method Editor Information Letter*.



## 2 UNICORN 6.x compatibility

### In this chapter

<b>Section</b>		<b>See page</b>
2.1	Supported operating systems	18
2.2	Supported SQL Server and databases	19
2.3	Instrument configurations compatibility	20

## 2.1 Supported operating systems

UNICORN versions	Windows Operating System		
	Win XP SP3	Win 7 SP1	OS Language
6.4.1 SP4	N/A	32/64-bit	US English
6.4.1 SP3	N/A	32/64-bit	US English
6.4.1 SP2	N/A	32/64-bit	US English
6.4.1 SP1	N/A	32/64-bit	US English
6.4.1	N/A	32/64-bit	US English
6.4.0	N/A	32/64-bit	US English
6.3.2 SP3	N/A	32/64-bit	US English
6.3.2 SP2	Yes	32/64-bit	US English
6.3.2 SP1	Yes	32/64-bit	US English
6.3.2	Yes	32/64-bit	US English
6.3.1	Yes	32/64-bit	US English
6.3	Yes	32/64-bit	US English
6.2	Yes	32/64-bit	US English

## 2.2 Supported SQL Server and databases

UNICORN version	SQL Server Express Edition	Database version	SQL Server tested combination				
			2008 R2	2012 SP1	2014 SP1	2016 SP2	2017
6.4.1 SP4	2008 R2	<b>1.6.7</b> and 1.7.4	X	X	X		
6.4.1 SP3	2008 R2	<b>1.6.7</b>	X	X	X		
6.4.1 SP2	2008 R2	<b>1.6.7</b>	X	X	X		
6.4.1 SP1	2008 R2	<b>1.6.7</b>	X				
6.4.1	2008 R2	<b>1.6.7</b>	X			X	X
6.4	2008 R2	<b>1.6.7</b>	X				
6.3.2 SP3	2008 R2	<b>1.6.6</b> , 1.6.7, and 1.7.4	X	X	X		
6.3.2 SP2	2008 R2	<b>1.6.6</b> and 1.6.7	X	X	X		
6.3.2 SP1	2008 R2	<b>1.6.6</b> and 1.6.7					
6.3.2	2008 R2	<b>1.6.6</b> and 1.6.7				X	X
6.3.1	2008 R2	<b>1.6.6</b>					
6.3	2008 R2	<b>1.6.5</b>					
6.2	2008 R2	<b>1.6.4</b>					

## 2.3 Instrument configurations compatibility

### In this section

<b>Section</b>	<b>See page</b>
2.3.1 ÄKTA avant 150	21
2.3.2 ÄKTA avant 25	22
2.3.3 ÄKTA pure 150	23
2.3.4 ÄKTA pure 25	24
2.3.5 Several ÄKTA systems, ReadyToProcess WAVE, and UniFlux	25
2.3.6 Legacy ÄKTA systems	26

## 2.3.1 ÄKTA avant 150

ÄKTA avant 150 IC	UNICORN version							
	6.1	6.2	6.3	6.3.1	6.3.2	6.3.2 AC	6.4	6.4.1
1.0.5.0	X	X	X	X	X	X	X	X
1.1.2.0	X	X	X	X	X	X	X	X
1.2.5.0	X	X	X	X	X	X	X	X
1.3.1.0	X	X	X	X	X	X	X	X
1.4.0.0	X	X	X	X	X	X	X	X
1.5.1.0	X	X	X	X	X	X	X	X
1.6.0.0	X	X	X	X	X	X	X	X
2.0.2.1				X	X	X	X	X
2.1.0.2				X <sup>1</sup>	X <sup>1</sup>	X	X	X
2.2.0.0				X <sup>1</sup>	X <sup>1</sup>	X	X	X
2.3.0.0				X <sup>1</sup>	X <sup>1</sup>	X	X	X
2.4.0.4				X <sup>1</sup>	X <sup>1</sup>	X	X	X
2.5.0.0				X <sup>1</sup>	X <sup>1</sup>	X	X	X
2.6.0.3				X <sup>1</sup>	X <sup>1</sup>	X	X	X
2.7.0.5				X <sup>1</sup>	X <sup>1</sup>	X	X	X

<sup>1</sup> This instrument configuration version does not support Win XP even though UNICORN 6.3.2 does.

## 2.3.2 ÄKTA avant 25

ÄKTA avant 25 IC	UNICORN version								
	6.0	6.1	6.2	6.3	6.3.1	6.3.2	6.3.2 AC	6.4	6.4.1
1.0.9.0	X								
1.1.2.0	X								
1.2.0.0	X								
1.3.0.0	X								
1.4.0.2	X								
2.0.5.0		X	X	X	X	X	X	X	X
2.1.2.0		X	X	X	X	X	X	X	X
2.2.5.0		X	X	X	X	X	X	X	X
2.3.1.0		X	X	X	X	X	X	X	X
2.4.1.0		X	X	X	X	X	X	X	X
2.5.0.0		X	X	X	X	X	X	X	X
3.0.2.1						X <sup>1</sup>	X <sup>1</sup>	X	X
3.1.0.3						X <sup>1</sup>	X <sup>1</sup>	X	X
3.2.0.0						X <sup>1</sup>	X <sup>1</sup>	X	X
3.3.0.0						X <sup>1</sup>	X <sup>1</sup>	X	X
3.4.0.15						X <sup>1</sup>	X <sup>1</sup>	X	X
3.5.0.0						X <sup>1</sup>	X <sup>1</sup>	X	X
3.6.0.15						X <sup>1</sup>	X <sup>1</sup>	X	X
3.7.0.4						X <sup>1</sup>	X <sup>1</sup>	X	X

<sup>1</sup> This instrument configuration version does not support Win XP even though UNICORN 6.3.2 does.

### 2.3.3 ÄKTA pure 150

ÄKTA pure 150 IC	UNICORN version					
	6.3	6.3.1	6.3.2	6.3.2 AC	6.4	6.4.1
1.0.0.1	X	X	X	X	X	X
1.2.0.3			X	X	X	X
1.3.0.1			X	X	X	X
1.4.1.2			X	X	X	X
1.5.0.0			X	X	X	X
1.6.0.2			X	X	X	X
1.7.0.6			X	X	X	X
1.8.0.3			X	X	X	X

## 2.3.4 ÄKTA pure 25

ÄKTA pure 150 IC	UNICORN version					
	6.3	6.3.1	6.3.2	6.3.2 AC	6.4	6.4.1
1.0.1.2	X	X	X	X	X	X
1.1.4.0	X	X	X	X	X	X
1.2.0.0	X	X	X	X	X	X
1.4.0.3			X	X	X	X
1.5.0.2			X	X	X	X
1.6.1.2			X <sup>1</sup>	X <sup>1</sup>	X	X
1.7.0.1			X <sup>1</sup>	X <sup>1</sup>	X	X
1.8.0.6			X <sup>1</sup>	X <sup>1</sup>	X	X
1.9.0.11			X <sup>1</sup>	X <sup>1</sup>	X	X
1.10.0.6			X <sup>1</sup>	X <sup>1</sup>	X	X
1.12.0.6			X <sup>1</sup>	X <sup>1</sup>	X	X

<sup>1</sup> This instrument configuration version does not support Win XP even though UNICORN 6.3.2 does.



## 2.3.5 Several ÄKTA systems, ReadyToProcess WAVE, and UniFlux

Instrument configuration	UNICORN version						
	6.2	6.3	6.3.1	6.3.2	6.3.2 AC	6.4	6.4.1
(ÄKTAprocess) 29165862_AA_W230(1.0.0.0)	X	X	X	X	X	X	X
Custom bioprocess				X	X	X	X
<b>(ReadyToProcess)</b> WAVE 25/Xuri 1.0.3.0					X	X	X
WAVE 25/Xuri 2.0.4.0					X	X	X
WAVE 25/Xuri 2.0.7.0						X	X
WAVE 25/Xuri 2.0.8.0						X	X
WAVE 25/Xuri 2.0.9.0						X	X
UniFlux				X	X	X	X

## 2.3.6 Legacy ÄKTA systems

Instrument configuration	UNICORN version						
	6.2	6.3	6.3.1	6.3.2	6.3.2 AC	6.4	6.4.1
(ÄKTApilot) PiloX_133_W253_1.0.0.2	X	X	X	X	X	X	X
<b>(ÄKTAexplorer)</b> E10_414_1.0.0.0						X	X
E10F414_1.0.0.0						X	X
E100_444_1.0.0.0						X	X
E100F444_1.0.0.0						X	X
<b>(ÄKTApurifier/Basic)</b> P10F414_1.0.0.0						X	X
P10_415_1.0.0.0						X	X
P100_415_1.0.0.0						X	X
P100F415_1.0.0.0						X	X
<b>(ÄKTApurifier UPC)</b> P10_UPC_414_1.0.0.0						X	X
P10F_UPC_414_1.0.0.0						X	X
P100_UPC_414_1.0.0.0						X	X
P100F_UPC_414_1.0.0.0						X	X

## 3 Support for RTU and RTU Win10 Image

UNICORN version											
	7.1 and earlier	7.2	7.3	7.4	7.5	7.6	7.7	7.8	7.9	7.10	7.11
RTU	N/A	X	X	X	X	X	X	X	X	X	X

### UNICORN, IC, and RTU Win10 Image combinations for upgrade

RTU Win10 Image	UNICORN version	IC	Valid combination?
v1.15 or earlier	7.6 or later	Signed by Cytiva	Yes
v1.15 or earlier	7.6 or later	Signed by GE	Yes
v1.15 or earlier	7.5 or earlier	Signed by Cytiva	No, upgrade to 7.6 or later
v1.15 or earlier	7.5 or earlier	Signed by GE	Yes
v1.15 or later	7.6 or later	Signed by Cytiva	Yes
v1.15 or later	7.6 or later	Signed by GE	Yes
v1.15 or later	7.5 or earlier	Signed by Cytiva	Yes
v1.15 or later	7.5 or earlier	Signed by GE	Yes

For Instrument Configurations with older Cytiva certificates, an update of the RTU Win10 Image is needed (for image versions 2.0 or earlier): RTU Win10 Image SP1 must be installed. For more information, see the specific IC version affected.

# 4 Modifications history

This chapter describes the history of modifications done on the current version of this document with respect to information released on previous versions.

<b>UNICORN version</b>	<b>Type of modification</b>	<b>Instrument Configuration</b>	<b>Modification date</b>
7.9, 7.10	Compatibility removed	ÄKTA avant 150 2.4.0.4, 2.5.0.0, 2.6.0.3, 2.7.0.5, 3.0.0.8, and 3.1.0.30	June 2024
7.9, 7.10	Compatibility removed	ÄKTA avant 25 3.4.0.15, 3.5.0.0, 3.6.0.15, 3.7.0.4, 4.0.0.8, and 4.1.0.30	June 2024
7.9, 7.10	Compatibility removed	ÄKTA pure 150 1.6.0.2, 1.7.0.6, 1.8.0.3, 2.0.0.4, 2.1.0.13, and 2.2.0.3	June 2024
7.9, 7.10	Compatibility removed	ÄKTA pure 25 1.8.0.6, 1.9.0.11, 1.10.0.6, 1.12.0.6, 1.13.0.63, 2.0.0.20, 2.1.0.1, 2.2.0.27, and 2.3.0.3	June 2024
7.9, 7.10	Compatibility removed	ÄKTA pilot 600 1.0.0.53, 1.1.0.6, and 1.2.0.2	June 2024
7.9, 7.10	Compatibility removed	ÄKTA go 1.0.0.17, 2.0.0.44, 2.1.0.0, and 2.2.0.1	June 2024

Page intentionally left blank

**Give feedback on this document**

Visit [cytiva.com/techdocfeedback](https://cytiva.com/techdocfeedback) or scan the QR code.



## cytiva.com/unicorn

Cytiva and the Drop logo are trademarks of Life Sciences IP Holdings Corporation or an affiliate doing business as Cytiva.

ÄKTA, ÄKTA avant, ÄKTA crossflow, ÄKTA go, ÄKTA oligosynt, ÄKTA pcc, ÄKTA pilot, ÄKTAprocess, ÄKTA process, ÄKTA pure, ÄKTA ready, ÄKTA readyflux, OligoPilot, OligoProcess, ReadyToProcess, ReadyToProcess WAVE, UNICORN, UniFlux, and Xuri are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Any other third-party trademarks are the property of their respective owners.

© 2020–2024 Cytiva

UNICORN © 2020–2024 Cytiva

Any use of software may be subject to one or more end user license agreements, a copy of, or notice of which, are available on request.

For local office contact information, visit [cytiva.com/contact](https://cytiva.com/contact)

29258394 BL V:23 12/2024