

UNICORN™

Computer Specifications

Minimum computer specifications

	UNICORN™ Client	Database Server	Workstation installation	E-License Server
Free disk space	16 GB	10 GB	27 GB	500 MB
Available RAM	4 GB	4 GB	8 GB	2 GB
Disc format	NTFS	NTFS	NTFS	NTFS
Architecture	Quad-core processor or 4 logical processors	Quad-core processor or 4 logical processors	Quad-core processor or 4 logical processors	Multi-core processor



IMPORTANT

Two network cards are necessary for the computers running the Instrument Server.

**Give feedback on this document**

Visit 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.

UNICORN is a trademark 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.

© 2020-2025 Cytiva

UNICORN © 2020-2025 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

29283583 AE V:4 01/2025