

Amersham ECL Rainbow Marker - Low Range

Product Specification Sheet

Introduction

Product code

RPN755E

About

$M_r = 3\,500 - 40\,000$

Important

Read these instructions carefully before using the products.

Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.

Storage

Store at -15°C to -30°C . Stable for at least 3 months when stored under recommended conditions.

Concentration

Approx. 1 mg/mL of protein.

Pack size

250 μL , sufficient for for 50 minigel loadings when used under recommended conditions.

Description

Amersham™ ECL™ Rainbow™ Marker - Low Range is a mixture of individually colored proteins of defined size from Cytiva. Purified proteins are combined to produce bands of equal color intensity and even spacing when separated on a polyacrylamide gel as described by Laemmli (1), Schagger and von Jagow (2), Swank and Munkres (3), Weber and Osborn (4).

Form

Supplied ready to use in 30% glycerol and sample buffer containing mercaptoethanesulphonic acid (MESNA) as reducing agent (5).

Molecular weight (Da)	Color
38 000	Blue
31 000	Orange
24 000	Green
17 000	Blue
12 000	Red

Molecular weight (Da)	Color
8500	Yellow
3500	Blue

Usage

Recommended minimum loadings are as follows:

8 × 10 cm gels: 5 μL of Amersham ECL Rainbow Marker - Low Range.

20 × 20 cm gels: 10 μL of Amersham ECL Rainbow Marker - Low Range.

Step	Action
1	Remove the marker from storage at -15°C to -30°C and allow to equilibrate to room temperature. A precipitate of SDS may form on storage at -15°C to -30°C . If necessary briefly warm the solution at 37°C to dissolve the precipitate.
2	Mix well and load the required volume of markers directly on to the gel.

More technical help, tips, and best practices can be found in the handbook *Western Blotting Principles and Methods* from Cytiva (Product code 28999897).

Typical result RPN755E

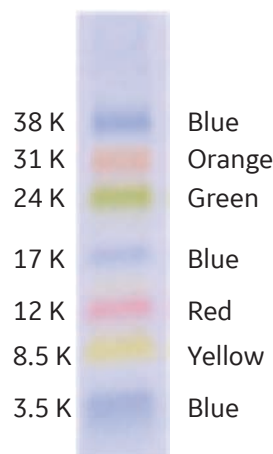


Fig 1. 4–20% Tris-Glycine gradient SDS-PAGE gel. 18% Tris-Glycine SDS-PAGE gel Electrophoresis was performed for 90 minutes at 125 V.

Quality control

Each batch of Amersham ECL Rainbow Marker - Low Range is assessed for color intensity and band integrity on an 18% Tris-Glycine SDS-PAGE mini-gel.

24 k Green band

In some gel/buffer systems the mobility of this band may differ from that quoted using a Tris/Glycine/SDS buffer.

Measurement of protein sizes

The sizes of the labeled proteins have been determined by interpolation from a standard curve of Rf values of known molecular weight recombinant proteins on a 4–20% Tris-Glycine gradient SDS-PAGE gel.

Related products

Amersham ECL DualVue™ Western Blotting RPN810

Markers (M_r = 15 000 – 150 000)

Amersham ECL Plex™ Fluorescent Rainbow RPN850E, RPN851E

Markers (M_r = 12 000 – 225 000)

Amersham ECL Rainbow Marker - High Range, (M _r = 12 000 – 225 000)	RPN756E (10 protein ladder)
---	--------------------------------

Amersham ECL Rainbow Marker - Full Range, (M _r = 12 000 – 225 000)	RPN800E (10 protein ladder)
---	--------------------------------

References

1. Laemmli, U.K., *Nature* **227**, 681 (1970).
2. Schagger, H. and von Jagow, G., *Anal. Biochem.* **166**, 368 (1987).
3. Swank, R.T. and Munkres, K.D., *Anal. Biochem.* **39**, 462 (1971).
4. Weber, K. and Osborn, M., *J. Biol. Chem.* **244**, 4406 (1969).
5. Singh, R., *Biotechniques*. **17**, 263 (1994).

cytiva.com

Cytiva and the Drop logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Amersham, ECL, ECL DualVue, ECL Plex and Rainbow are trademarks of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

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

© 2020–2021 Cytiva

All goods and services are sold subject to the terms and conditions of sale of the supplying company operating within the Cytiva business. A copy of those terms and conditions is available on request. Contact your local Cytiva representative for the most current information.

For local office contact information, visit cytiva.com/contact

28925388 AE V:4 02/2021

