One-Step Lumitein™ Protein Gel Stain

A safe, ready-to-use luminescent protein gel stain. Because it requires no fixation or wash steps, it is a big improvement on the original LumiteinTM protein gel stain.



Product attributes

Call us: 800-304-5357

Excitation/Emission

280, 460 (broad)/610 nm

Email: btinfo@biotium.com

Product Description

One-Step Lumitein[™] Protein Gel Stain is a ready-to-use luminescent red protein gel stain. It is a dramatically improved version of our original Lumitein[™] protein gel stain in terms of convenience and safety. One-Step Lumitein[™] can detect as little as approximately 1-10 ng of protein per band.

- One-step staining procedure, no fixation required
- Non-toxic formulation is safe for easy disposal
- Sensitive protein detection down to 1-10 ng
- Image with a UV transilluminator, a blue light transilluminator, or laser-based gel scanner
- 1 L and 4 L sizes
- Store at 4°C

Fast and simple procedure

One-Step LumiteinTM gel staining requires only a single 5-60 minute staining step without fixation. Destaining is optional. One-Step LumiteinTM can detect as little as approximately 1-10 ng of protein per band depending on the staining method used, although staining intensity varies between proteins. Staining is fully compatible with mass spectrometry and Edman-based sequencing. Staining can be imaged using a UV transilluminator with EtBr or SYBR filter set, a blue light transilluminator, or a laser-based gel scanner.

One-Step Lumitein™ is non-toxic

One-Step LumiteinTM Protein Gel Stain offers safer handling and ease of disposal compared to other protein gel stains, because it is an aqueous-based solution that does not contain hazardous methanol or acetic acid. One-Step stains are certified under CCR Title 22 as non-toxic to the environment for drain disposal after a simple pH neutralization step. Furthermore, One-Step LumiteinTM costs significantly less than other fluorescent protein gel stains. One-Step LumiteinTM Protein Gel Stain is available in 1 liter bottle or 4 liter Cubitainer® sizes.

Please visit our <u>Protein Gel Stains technology page</u> to learn about our other safe and sensitive protein gel stains such as <u>One-Step LumiteinTM UV</u>, which is optimized for use with a UV transilluminator, and <u>One-Step BlueTM</u>, a rapid, easy-to-use, non-toxic alternative to Coomassie staining for visible blue staining.

Product	Replacement for	Catalog number	Unit size	Product protocol	Safety report
One-Step Blue®	Coomassie Brilliant Blue stains	21003-1L	1L	<u>PI-21003</u>	One-Step Gel Stains Safety Report
One-Step Lumitein™	SYPRO® Ruby Protein Gel Stain	21004-1L	1L	<u>PI-21004</u>	
		21004-4L	4L		
One-Step Lumitein™ UV	Oriole™ and Flamingo™ Protein Gel Stains	21005-1L	1L	<u>PI-21005</u>	
		21005-4L	4L		

One-Step Blue is a registered trademark of Biotium. Lumitein and its related technologies are covered by US and international patents. SYPRO is a registered trademark of Molecular Probes, Inc. Oriole and Flamingo are trademarks of Bio-Rad Laboratories. Cubitainer is a registered trademark of Hedwin Corporation.

This datasheet was generated on November 9, 2025 at 09:03:58 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/one-step-lumitein-protein-gel-stain/