



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## 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 Lumitein™ protein gel stain.



### Product attributes

Excitation/Emission	280, 460 (broad)/610 nm
---------------------	-------------------------

## 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 Lumitein™ gel staining requires only a single 5-60 minute staining step without fixation. Destaining is optional. One-Step Lumitein™ 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 Lumitein™ 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 Lumitein™ costs significantly less than other fluorescent protein gel stains. One-Step Lumitein™ Protein Gel Stain is available in 1 liter bottle or 4 liter Cubitainer® sizes.

Please visit our [Protein Gel Stains technology page](#) to learn about our other safe and sensitive protein gel stains such as [One-Step Lumitein™ UV](#), which is optimized for use with a UV transilluminator, and [One-Step Blue™](#), 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
<a href="#">One-Step Blue®</a>	Coomassie Brilliant Blue stains	21003-1L	1L	<a href="#">PI-21003</a>	<a href="#">One-Step Gel Stains Safety Report</a>
<a href="#">One-Step Lumitein™</a>	SYPRO® Ruby Protein Gel Stain	21004-1L	1L	<a href="#">PI-21004</a>	
21004-4L	4L				
<a href="#">One-Step Lumitein™ UV</a>	Oriole™ and Flamingo™ Protein Gel Stains	21005-1L	1L	<a href="#">PI-21005</a>	
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 January 9, 2026 at 04:43:02 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/one-step-lumitein-protein-gel-stain/>