



Glowing products for science

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

# One-Step Lumitein™ UV Protein Gel Stain

A ready-to-use luminescent protein gel stain optimized for UV-based fluorescence imaging systems. Non-toxic, with no fixation required.



## Product attributes

Excitation/Emission	288/603 nm
---------------------	------------

## Product Description

One-Step Lumitein™ UV Protein Gel Stain is a ready-to-use luminescent red protein gel stain optimized for use on UV transilluminators. With no fixation or wash steps required, One-Step Lumitein™ UV is a great choice for convenience and safety. One-Step Lumitein™ UV 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
- Optimized for use with a UV transilluminator
- 1 L and 4 L sizes
- Store at 4°C

## Fast and simple procedure

One-Step Lumitein™ UV gel staining requires only a single 5-30 minute staining step without fixation. Destaining is optional. One-Step Lumitein™ UV 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. Gels stained with One-Step Lumitein™ UV Protein Gel Stain can be visualized with a variety of different UV-based fluorescence imaging systems.

The results obtained with One-Step Lumitein™ UV are comparable to those with Oriole™ Fluorescent Gel Stain; however, One-Step Lumitein™ UV has a more convenient protocol, does not contain hazardous solvents, and does not cause gel shrinkage. Furthermore, One-Step Lumitein™ costs significantly less than other fluorescent protein gel stains. One-Step Lumitein™ UV Protein Gel Stain is available in 1 liter bottle or 4 liter Cubitainer® sizes.

## One-Step Lumitein™ UV is non-toxic

One-Step Lumitein™ UV Protein Gel Stain offers safer handling and ease of disposal, 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.

Please visit our [Protein Gel Stains technology page](#) to learn about our other safe and sensitive protein gel stains such as [One-Step Lumitein™](#) 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.

## References

Nature Biotechnology (2018), doi:10.1038/nbt.4173

This datasheet was generated on December 21, 2025 at 09:07:09 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/one-step-lumitein-uv-protein-gel-stain/>