# One-Step Lumitein™ UV Protein Gel Stain

TO THE STATE OF TH

Product attributes

Excitation/Emission

Call us: 800-304-5357 Email: btinfo@biotium.com

288/603 nm

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

## **Product Description**

One-Step Lumitein<sup>TM</sup> 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<sup>TM</sup> UV is a great choice for convenience and safety. One-Step Lumitein<sup>TM</sup> 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<sup>TM</sup> UV gel staining requires only a single 5-30 minute staining step without fixation. Destaining is optional. One-Step Lumitein<sup>TM</sup> 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<sup>TM</sup> UV Protein Gel Stain can be visualized with a variety of different UV-based fluorescence imaging systems.

The results obtained with One-Step Lumitein<sup>TM</sup> UV are comparable to those with Oriole<sup>TM</sup> Fluorescent Gel Stain; however, One-Step Lumitein<sup>TM</sup> UV has a more convenient protocol, does not contain hazardous solvents, and does not cause gel shrinkage. Furthermore, One-Step Lumitein<sup>TM</sup> costs significantly less than other fluorescent protein gel stains. One-Step Lumitein<sup>TM</sup> 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 <a href="CCR Title 22">CCR Title 22</a> as non-toxic to the environment for drain disposal after a simple pH neutralization step.

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 Lumitein<sup>TM</sup></u> and <u>One-Step Blue<sup>TM</sup></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	PI-21003	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.

### References

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

This datasheet was generated on November 6, 2025 at 07:05:10 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/one-step-lumitein-uv-protein-gel-stain/