



Glowing products for science

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

4X Protein Loading Buffer with Orange Tracking Dye

A convenient 4X sample loading buffer for protein gel electrophoresis. The loading buffer is optimized for use with denaturing SDS-PAGE.



Product attributes

Product Description

4X Protein Loading Buffer with Orange Tracking Dye is a convenient loading buffer for protein gel electrophoresis. The loading buffer contains an orange tracking dye that is recommended for fluorescent gel and western blot applications because it avoids unwanted background fluorescence caused by blue tracking dyes in far-red and near-infrared detection channels. The loading buffer is formulated for denaturing SDS-PAGE and is compatible with both visible fluorescent and near-IR fluorescent gel and western blot imaging.

Learn more about Biotium's other products for [western blotting](#) including reversible [VersaBlot™ Total Protein Normalization Kits](#) and [Peacock™ Prestained Protein Markers](#).

This datasheet was generated on June 23, 2026 at 05:11:10 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/4x-protein-loading-buffer-with-orange-tracking-dye/>