



Glowing products for science

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

AccuOrange™ Protein Quantitation Kit

Highly sensitive, fluorescence-based quantitation of purified proteins.



Product attributes

Excitation/Emission	480/598 nm
Storage Conditions	Store at room temperature, Do not refrigerate, Protect dye component from light

Product Description

AccuOrange™ Protein Quantitation Kit is a highly sensitive fluorescence-based assay for quantifying purified protein samples in 96-well format.

- Highly sensitive: detect 0.1-15 ug/mL protein
- Excellent linearity and low variability
- Stable fluorescence signal
- Compatible with reducing agents and other additives
- For use with purified protein or antibody samples
- Ex/Em: 480/598 nm

The detection range of the assay is 0.1-15 ug/mL protein. AccuOrange™ is much more sensitive than traditional protein quantitation assays such as BCA, Bradford and Lowry, and shows superior linearity and reproducibility compared to the NanoOrange® protein quantitation assay. The assay shows minimal variability between different proteins, and has stable fluorescence signal for up to 16 hours.

AccuOrange™ is recommended for quantifying purified protein or antibody samples. The tolerance of the AccuOrange™ assay to salts, buffers, detergents, and other chemicals is listed in the product protocol. The AccuOrange™ assay has low tolerance for non-ionic detergents, and is not recommended for use with cell lysates containing Triton™ X-100, sodium deoxycholate, CHAPs, or other non-ionic detergents. The assay can tolerate up to 0.01% SDS (final concentration in assay).

NanoOrange is a registered trademark of Thermo Fisher Scientific. Triton is a trademark of The Dow Chemical Company.

This datasheet was generated on June 19, 2026 at 08:42:17 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/accuorange-protein-quantitation-kit/>