



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).

Comparison of AccuOrange™ with other protein quantitation assays

Assay	Detection Range	Comments
AccuOrange™	0.1-15 ug/mL	Fluorescence-based (480/598 nm) Highly linear Signal stable for at least 16 hours Compatible with reducing agents Not compatible with detergents
NanoOrange®	0.1-10 ug/mL	Fluorescence-based (470/570 nm) Non-linear Fluorescence stable for 6 hours Compatible with reducing agents Not compatible with detergents
Modified Lowry	1-1500 ug/mL	Absorbance-based (750 nm) Non-linear Not compatible with reducing agents or detergents
BCA	20-2000 ug/mL	Absorbance-based (562 nm) Highly linear Signal not stable over time Not compatible with reducing agents Compatible with detergents
Bradford (Coomassie®)	50-500 ug/mL	Absorbance-based (595 nm) Signal not stable over time Non-linear Compatible with reducing agents Not compatible with detergents
Pierce® 660 nm	50-2000 ug/mL	Absorbance-based (660 nm) Non-linear Compatible with reducing agents & detergents
A ₂₈₀	50-2000 ug/mL	Absorbance-based (280 nm) High protein-protein variability Contaminants and detergents can affect results

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

This datasheet was generated on January 14, 2026 at 06:16:56 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/accuorange-protein-quantitation-kit/>