

Product Information

AccuOrange™ Protein Quantitation Kit

Catalog Number: 30071-T (200 assays), 30071 (2000 assays)

Kit Contents

Component	30071-T 200 assays	30071 2000 assays
AccuOrange buffer, 10X (contains 2 mM sodium azide)	5 mL 30071A-T	50 mL 30071A
AccuOrange dye, 500X in DMSO	100 uL 30071B-T	1 mL 30071B
Bovine serum albumin (BSA) standard, 2 mg/mL (contains 2 mM sodium azide)	100 uL 99990-T	1 mL 99990

Number of assays based on 96-well format.

Storage and Handling

Store at room temperature. Do not refrigerate AccuOrange buffer. Protect AccuOrange dye from light. Product is stable for at least 1 year from date of receipt when stored as recommended. If precipitate forms in Assay Buffer, heat to 37°C and swirl to re-dissolve. No data are available on the safety of AccuOrange dye. Handle using universal laboratory safety precautions and dispose of the dye as chemical waste.

Spectral Properties

Ex/Em 480/598 nm with BSA in 1X assay buffer (Fig. 1)

Product Description

AccuOrange™ Protein Quantitation Kit is a highly sensitive fluorescence-based assay for quantitating purified protein samples in 96-well format. 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 (Figure 2). The assay shows minimal variability between different proteins. After protein samples are heated with the dye, the fluorescence signal is stable for up to 16 hours at room temperature.

AccuOrange is recommended for quantitating purified protein or antibody samples. The tolerance of the AccuOrange assay to salts, buffers, detergents, and other chemicals is shown in Table 1. 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).

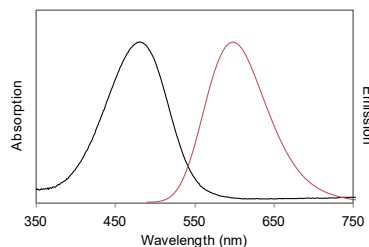


Figure 1. Absorbance and emission spectra of AccuOrange dye with BSA in 1X AccuOrange assay buffer.

Assay Protocol

1. Prepare 1X AccuOrange buffer by diluting 10X buffer (component 30071A) 1:10 in dH₂O. For example, add 1 mL of 10X assay buffer to 9 mL dH₂O. 1X assay buffer can be stored at room temperature.
2. Immediately before use, prepare AccuOrange working solution by diluting 500X AccuOrange dye (component 30071B) 1:500 in 1X AccuOrange buffer. For example, add 10 uL of 500X AccuOrange dye to 5 mL of 1X assay buffer.

Note: You will need about 3 mL working solution for each standard curve (see Table 1) and 250 uL working solution for each sample.
3. Prepare unknown samples by adding up to 10 uL of sample to 250 uL AccuOrange working solution.

Note: You may wish to perform a few different dilutions of your unknown sample. See Table 2 for the assay tolerance levels for buffer components and potential contaminants. Sample dilution may reduce the concentration of interfering substances to tolerable levels.

4. Prepare a protein standard curve by performing serial dilutions of the BSA standard in AccuOrange working solution as shown in Table 1.

Note: the standards will be measured at the concentrations shown in Table 1, without further dilution.

5. Heat samples and standards to 90°C-95°C for 10 minutes, protected from light. Samples can be heated in microcentrifuge tubes in a water bath or heat block. Screw cap tubes or cap locks are recommended to prevent caps from popping open during heating.

Note: Alternatively, samples can be heated in a multiwell plate with a heat-resistant seal using a thermocycler with a heated lid. Scale all volumes in the assay proportionally to fit the well volume of the plate. Centrifuge the plate to collect any condensation that forms on plate seal after step 6.

6. Remove samples from heat and allow to cool to room temperature for 20 minutes, protected from light. Centrifuge tubes briefly to collect any condensation from caps and vortex to mix.

7. Transfer 200 uL of each standard or sample to 96-well microplate wells to read on a fluorescence microplate reader. Measure fluorescence with excitation/emission at 480/598 nm.

Note: Alternatively, samples can be transferred to a fluorescence cuvette and measured using a spectrofluorometer. If more than 200 uL is required for measurement, scale all volumes in the assay proportionally.

Table 1. Preparation of BSA standards.

	Volume of BSA solution	Volume of working solution	Final BSA concentration (ug/mL)
A	7.5 uL BSA std (2 mg/mL)	1 mL	15
B	333 uL solution A	167 uL	10
C	250 uL solution B	250 uL	5
D	250 uL solution C	250 uL	2.5
E	200 uL solution D	300 uL	1
F	250 uL solution E	250 uL	0.5
G	250 uL solution F	250 uL	0.25
H	100 uL solution G	150 uL	0.1
I	0 mL	250 uL	0

Table 2. Assay tolerance levels for interfering substances

Compound	Maximum tolerable concentration (final concentration in assay)
SDS	0.01%
Triton X-100	Below 0.001%
Tween 20	Below 0.001%
CHAPS	Below 0.001%
Sodium deoxycholate (DOC)	Below 0.001%
Urea	10 mM
DTT	100 mM
β ME	0.1%
EDTA	1 mM
Sucrose	10 mM (0.34%)
Glycerol	1%
PBS	0.02X
NaCl	1 mM
CaCl_2	0.01 mM
MgCl_2	0.2 mM
ZnCl_2	0.2 mM
Ammonium sulfate	1 mM
Sodium azide	2 mM
Imidazole	50 mM
Tris	1 mM
Glycine	100 mM
Amino acids	10 $\mu\text{g/mL}$
DNA	10 $\mu\text{g/mL}$

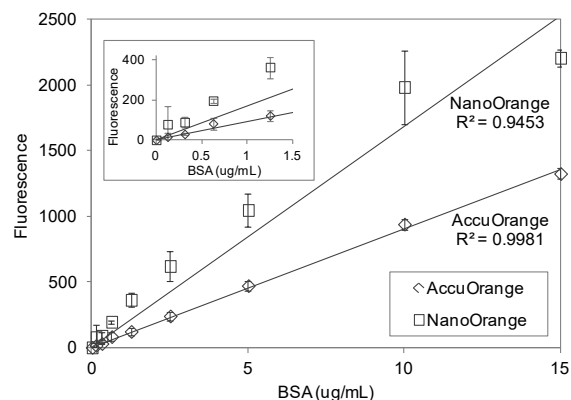


Figure 2. BSA titration assayed using AccuOrange Protein Quantitation Kit or NanoOrange Protein Quantitation Kit from Life Technologies according to manufacturer's protocol and read on a microplate reader at the recommended wavelengths for each assay. Inset shows the lower end of the curve. Error bars represent standard deviation of the mean for triplicate samples.

Related Products

Catalog number	Product
21003-1L	One-Step Blue™ Protein Gel Stain
21004	One-Step Lumitein™ Protein Gel Stain
21005-1L	One-Step Lumitein™ UV Protein Gel Stain
31060	AccuBlue™ NextGen dsDNA Quantitation Kit
31028	AccuClear™ Ultra High Sensitivity dsDNA Quantitation Kit with 7 DNA Standards
31006	AccuBlue™ High Sensitivity dsDNA Quantitation Kit
31007	AccuBlue™ Broad Range dsDNA Quantitation Kit
31066	AccuGreen™ High Sensitivity dsDNA Quantitation Kit (for Qubit)
41003	GelRed™ Nucleic Acid Gel Stain, 10,000X in water
31041-T	Forget-Me-Not™ qPCR Master Mix, trial size
31043-T	Forget-Me-Not™ Universal Probe Master Mix, trial size

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF™ dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

NanoOrange is a registered trademark of Thermo Fisher.

AccuOrange is a trademark of Biotium, Inc.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.