Ubiquitin (human), (fluorescein labeled)

The covalent attachment of ubiquitin to proteins (ubiquitinylation) and their subsequent proteasomal degradation plays a fundamental role in the regulation of cellular function through biological events involving cell cycle, differentiation, immune responses, DNA repair, chromatin structure, and apoptosis.

Citations: 4

View Online »

Ordering Information

Order Online »

BML-UW1240-0100

100µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Stable for at least

1 year after receipt when stored at -20°C. Reagent is light sensitive; store in the dark or in amber vials. Stable to freeze/thaw but this should be minimized wherever possible.

Handling Protect from light.

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Notes For use in ubiquitinylation studies.

Formulation Lyophilized.

MW ~9.5kDa

Purity ≥95% (SDS-PAGE)

Purity Detail Purified by multi-step chromatography.

Reconstitution Reconstitute in minimum volume of DMSO (5-10µL)

followed by dilution to required concentration with aqueous

buffer of choice (e.g. 1mg/mL in PBS).

Source Produced in *E. coli*.

Technical Info / Product Notes

Fluorescein labelled ubiquitin is a useful reagent for studying ubiquitinylation *in vitro*, where it represents a readily detectable alternative to radio-iodinated ubiquitin. Fluorescein labelled ubiquitin is activated by ubiquitin E1, is a competent substrate for known ubiquitin E2/E3 enzymes (thioester bond formation) and can be conjugated to ubiquitin target proteins (isopeptide bond formation).

This product contains a number of multiply fluoresceinlabelled ubiquitin species, with [fluorescein]₂₋₄ubiquitin being the predominant species.

Typical thioester assay set-up: 2.5μM Fluorescein-Ub (BML-UW1240), 100nM Ub-E1 (BML-UW9410), 2.5μM E2 (e.g. His₆-UbcH5a, BML-UW9050), 1U inorganic pyrophosphatase, 5mM Mg-ATP in 20mM TRIS-HCl, pH 7.5, 1mM DTT. Incubate at 37°C for 30-60 mins. Analyse by SDS-PAGE followed by fluorescence detection.

P0CG47 (UBB), P0CG48 (UBC), P62979 (RPS27A), P62987 (UBA52)

UniProt ID

Last modified: May 29, 2024

uk@enzolifesciences.com