

Alkyne-PEG4-FLAG

Alkyne-PEG4-FLAG enables the FLAG-tag (DYKDDDDK) attachment to any Azide-functionalized molecule via Cu(I)-catalyzed terminal Alkyne-Azide Click Chemistry (CuAAC). The resulting FLAG-tagged molecules can subsequently be detected by an anti-FLAG antibody that is either immobilized onto a matrix (for purification) or coupled with a fluorescent dye or reporter enzyme for direct or indirect detection, respectively. Molecule solubility and efficient FLAG-tag detection is ensured by the integrated PEG-linker.

Ordering Information

[Order Online »](#)

JBS-CLK-088L	5x0.5µmol
--------------	-----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance white to off-white solid

Formula $C_{62}H_{91}N_{13}O_{28}S$

MW 1498.53

Purity $\geq 95\%$ (HPLC)

Solubility water, PBS (up to 30 mM tested)

Technical Info / Product Notes For the Original Manufacturer's data sheet please [click here](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com