

FGFR substrate (biotinylated)

This peptide, Biotin-Ahx-KEAPEDLYKDFLTL, is derived from the KDR sequence and is a substrate for FGFR. The biotin allows peptide to be used in kinase assays with streptavidin-bound membranes.

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Fibroblast growth factor receptor substrate (biotinylated)
MW	2021.3
Purity	≥95%
Sequence	Biotin-Ahx-Lys-Glu-Ala-Pro-Glu-Asp-Leu-Tyr-Lys-Asp-Phe-Leu-Thr-Leu

Last modified: May 29, 2024



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