

I κ B kinase substrate (biotinylated)

This peptide sequence, RHDSGLDSMKD, is derived from the area surrounding serine 32 and serine 36 of I κ B α , which is phosphorylated by I κ B Kinase (IKK). The biotin allows the peptide to be used in IKK kinase assays with streptavidin-bound membranes.

Ordering Information

[Order Online »](#)

BML-P148-0001	1mg
---------------	-----

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

MW	1486.9
Sequence	Biotin-Arg-His-Asp-Ser-Gly-Leu-Asp-Ser-Met-Lys-Asp



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