

Axltide (biotinylated)

Protein kinase substrate

This peptide, biotin-Ahx-KKSRGDYMTMQIG, is derived from mouse insulin receptor 1 and is a substrate for Axl, DDR2, Mst1, and JAK2 kinases. 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

| | |
|------------|----------------------------------------------------------------|
| Appearance | Lyophilized solid. |
| MW | 1854.2 |
| Purity | ≥95% |
| Sequence | biotin-Ahx-Lys-Lys-Ser-Arg-Gly-Asp-Tyr-Met-Thr-Met-Gln-Ile-Gly |
| Solubility | Soluble in water. |

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