LRRKtide (biotinylated)

LRRK substrate

This peptide, Biotin-Ahx-RLGRDKYKTLRQIRQ, is derived from ezrin/moesin/radixin and is a substrate for LRRK kinase. The biotin allows the peptide to be used in kinase assays with streptavidin-bound membranes.

Ordering Information

Order Online »

BML-P268-0001 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

Appearance White to off-white powder.

Formulation Lyophilized.

Identity Determined by MS.

MW 2270.9

Purity ≥95% (HPLC)

Sequence Biotin-Ahx-Arg-Leu-Gly-Arg-Asp-Lys-Tyr-Lys-Thr-Leu-Arg-

Gln-Ile-Arg-Gln-OH

Solubility Soluble in water.

Synthetic. Source

Last modified: May 29, 2024



info-

eu@enzolifesciences.com