MAPKAPK2 substrate

This peptide, KKLNRTLSVA is derived from the N-terminus of glycogen synthase and is known to be a good substrate for MAPKAPK2.

Ordering Information

Order Online »

BML-P196-0001 1r

1mg

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended. Store, as supplied,

at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name MAP kinase-activated protein kinase 2 substrate, Lys-Lys-

Leu-Asn-Arg-Thr-Leu-Ser-Val-Ala

Appearance White to off-white solid.

Formulation Lyophilized.

Identity Determined by MS.

MW 1130.5

Purity ≥95% (HPLC)

Sequence Lys-Lys-Leu-Asn-Arg-Thr-Leu-Ser-Val-Ala

Soluble in water.

Source Synthetic.

Last modified: May 29, 2024



eu@enzolifesciences.com