Src substrate II (biotinylated)

This peptide, Biotin-Ahx-YIYGSFK, was derived through a random combinatorial peptide library method and is a specific and efficient substrate for the scr family of kinases. The $\rm K_m$ of YIYGSFK for p60c-src is 55 μ M.

Ordering Information

Order Online »

BML-P260-0001 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

Formulation Lyophilized.

MW 1217.0

Purity ≥95% (HPLC)

Sequence Biotin-Ahx-Tyr-Ile-Tyr-Gly-Ser-Phe-Lys-OH

Solubility Soluble in water.

Last modified: May 29, 2024



eu@enzolifesciences.com