Src substrate II

This peptide, YIYGSFK, was derived through a random combinatorial peptide library method and is a specific and efficient substrate for the scr family of kinases. The K $_{\rm m}$ of YIYGSFK for p60c-src is 55 μM .

Ordering Information

Order Online »

BML-P259-0001 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 877.3

Purity ≥95% (HPLC)

H-Tyr-Ile-Tyr-Gly-Ser-Phe-Lys-OH Sequence

Solubility Soluble in water.

Last modified: May 29, 2024

