# Peptide A

#### Src kinase inhibitor

Sequence is pp60<sup>V-src</sup> (137-157), which is located in the SH3/SH2 domain region. Non-competitive inhibitor of Src and EGFR. Selectively inhibits protein tyrosine phosphorylation in cell extracts. Peptide is not phosphorylated by Src. IC $_{50}$ =7.5  $\mu$ M for v-Src. Inhibitory target seems to be a catalytic negative regulatory domain of the tyrosine kinase.

## **Ordering Information**

Order Online »

**BML-P306-0500** 0.5mg

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**

**Alternative Name** pp60v-src (137-157)

MW 2482.8

≥97% **Purity** 

Val-Ala-Pro-Ser-Asp-Ser-Ile-Gln-Ala-Glu-Glu-Trp-Tyr-Phe-Sequence

Gly-Lys-Ile-Thr-Arg-Arg-Glu

Last modified: May 29, 2024



info-

eu@enzolifesciences.com