

# Ras inhibitory peptide

The peptide sequence corresponds to amino acid residues 1149-1158 of the guanine nucleotide-releasing factor hSos1, which is essential for the control of Ras activity. Blocks the binding of hSos1 to human Grb2, a protein that binds to activated receptor tyrosine kinases.

Citations: 1

[View Online »](#)

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** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** SOS-1 (human) (1149-1158)

**Appearance** White lyophilized powder.

**Formula**  $C_{53}H_{91}N_{19}O_{11}$

**MW** 1170.4

**Purity** ≥97%

**Sequence** H-Val-Pro-Pro-Pro-Val-Pro-Pro-Arg-Arg-Arg-OH

**Solubility** Soluble in water.

**Last modified: May 29, 2024**



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)