

HSP27tide

Synthetic peptide substrate derived from the mouse HSP27

Catalog # H31-58

Lot # K151-3

Product Description

The HSP27tide peptide sequence (RRLNRQLSVA-amide) is based on the mouse HSP27 (amino acid 80-85).

Molecular Weight

The theoretical molecular weight is 1211.4

Purity

The purity was determined to be 98.6% by HPLC analysis.

Formulation

1mg of peptide supplied as a lyophilized powder.

Storage and Stability

Store product at -20°C. For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

Reconstitution Protocol

Dilute peptide in distilled H₂O to a final concentration of 1mg/ml.

Related Kinases

HSP27tide can be utilized as a substrate for the following active protein kinases.

Product Name	Catalog Number
MAPKAPK2, Active	M40-11G
MAPKAPK3, Active	M41-10G

HSP27tide

Synthetic peptide substrate derived from the mouse HSP27

Catalog Number H31-58

Quantity 1 mg

Specific Lot Number K151-3

Purity 98.6%

Format 1 mg lyophilized powder

Stability 1yr At -20°C from date of shipment

Storage & Shipping Store product at -20°C. For optimal storage, aliquot diluted product into smaller quantities and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped at ambient temperature.