

ULKtide

Synthetic peptide evaluated as a substrate for ULK family

Catalog # U01-58

Lot # U1690-5

Product Description

The ULKtide synthetic peptide [YANWLAASIYLDGKKK] is routinely evaluated as a substrate for ULK family kinase assays.

Molecular Weight

The theoretical molecular weight is 1841.16.

Purity

The purity was determined to be 95.45% 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 20mM Tris-HCl, pH 7.5 solution to a final concentration of 1mg/ml.

ULKtide

Synthetic peptide evaluated as a substrate for ULK family

Catalog #	U01-58
Quantity	1 mg
Lot #	U1690-5
Purity	95.45%
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.