

PRODUCT DATA SHEET



PEPTIDE NAME: Calpain Inhibitor I (ALLN)

SEQUENCE: N-acetyl-Leu-Leu-Norleucinal
(N-acetyl-Leu-Leu-Nle-CHO)

SEQUENCE DESCRIPTION: Inhibitor of neutral cysteine proteases.

CATALOG NO.: P-120

LOT NO.: temp

MOLECULAR WEIGHT: 383.5

PURITY: >97% by TLC analysis.

AMINO ACID ANALYSIS AND IDENTITY: Confirms.

PHYSICAL APPEARANCE: Lyophilized solid.

SOLUBILITY: DMSO 25mg/ml

STORAGE: -20°C

OTHER: See attached Guide to Handling and Storing Peptides.

LITERATURE REFERENCES:

1. S.W. Sherwood *et al. Proc. Natl. Acad. Sci. USA* 1993 **90** 3353
2. X. Wang *et al. Cell* 1994 **77** 53

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.