

Product Data Sheet

Product Name: c-Myc Peptide Epitope

Catalog Number: AS-24194 (1 mg) Lot Number: See label on vial

AS-20590 (5 mg)

Sequence: H-Glu-Gln-Lys-Leu-lle-Ser-Glu-Glu-Asp-Leu-OH (3-letter code)

EQKLISEEDL (1-letter code)

Molecular Weight: 1203.3

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Initially try to dissolve a small amount of c-Myc peptide in water. If peptide does not dissolve, c-Myc peptide is freely soluble in 1% NH₄OH

Storage: c-Myc peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Reconstituted peptide can be aliquoted and stored at -20 °C or lower.

Description: c-Myc, the product of the c-myc proto-oncogene, is a helix-loop-helix leucine zipper phosphoprotein that regulates gene transcription in cell proliferation, cell differentiation and apoptosis. This peptide is a human c-myc epitope. Ref: Johnson, G. and L. Hersh, *J. Biol. Chem.* **269**, 7682 (1994); Chou, T. et al. *J. Biol. Chem.* **270**, 18961 (1995); Kato, G. and C. Dang, *FASEB J.* **6**, 3065 (1992). See also TNF-FRET 1 (#25349), TNF-FRET 2 (#27126).

For Research Use Only