



## Product Data Sheet

---

<b>Product Name:</b>	c-Myc Peptide Epitope	
<b>Catalog Number:</b>	AS-24194 (1 mg) AS-20590 (5 mg)	Lot Number: See label on vial
<b>Sequence:</b>	H-Glu-Gln-Lys-Leu-Ile-Ser-Glu-Glu-Asp-Leu-OH (3-letter code) EQKLISEEDL (1-letter code)	
<b>Molecular Weight:</b>	1203.3	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	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% NH<sub>4</sub>OH

**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*