

Product Data Sheet

Product Name: Caspase 1 (ICE) Substrate 3f, fluorogenic

Catalog Number: AS-25285-5 (5mg) Lot Number: See label on vial

Sequence: Ac - Trp-Glu-His-Cys - AFC (3-letter code)

Ac - WEHD - AFC (1-letter code)

Molecular Weight: 838.6

Peptide Purity: >95%

Appearance: Lyophilized yellow powder

Peptide Reconstitution: Initially try to dissolve a small amount of Caspase 1 (ICE) Substrate in water. If the peptide does not dissolve, use 1% NH OH.

Storage: Caspase 1 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: AFC (7-amino-4-trifluoromethylcoumarin)-derived caspase substrates are widely used for the fluorimetric detection of various caspase activities. Cleavage of AFC peptides by caspases generate strongly fluorescent AFC that is monitored fluorimetrically at 500-510 nm with excitation of 370-390 nm.

Published Citations:

Gurtu V, et al. *Anal Biochem* **251**, 98-102 (1997). Martins LM, et al. *J Biol Chem* **272**, 7421-30 (1997). Deveraux QL, et al. *Nature* **388**, 300-4 (1997). Xiang J, et al. *Proc Natl Acad Sci USA* **93**, 14559-63 (1996).

For Research Use Only