

**TEV Protease**Tobacco Etch Virus Protease, recombinant, *E. coli*

Cat. No.	Amount
LE-001	1.000 Units

E-N-L-Y-F-Q  **(S/G/A/M/C/H)**

TEV protease recognition site

Unit Definition: One unit of TEV Protease will cleave 3 µg of control substrate in a total reaction volume of 10 µl in 1 hour at 30 °C in 50 mM Tris-HCl (pH 7.5) with 0.5 mM EDTA and 1 mM DTT.

For general laboratory use.

Shipping: shipped on dry ice

Storage Conditions: store at -80 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 30 kDa

CAS#: 139946-51-3

Purity: >85 % (SDS-PAGE)

Form: liquid

Concentration: 10 Units/µl

Description:

TEV Protease is a highly sequence-specific cysteine protease that is found in the Tobacco Etch Virus (TEV). The TEV Protease recognition sequence is Glu-Asn-Leu-Tyr-Phe-Gln-(Ser) [ENLYFQ(S)] and cleavage occurs between the Gln and Ser residues. The amino acid in the P1' position can also be G, A, M, C, or H. TEV Protease is used to cleave affinity tags from fusion proteins. The protease has a His-tag for easy removal from a reaction using nickel affinity resins (#AC-501).

Content:**TEV Protease**

10 units/µl TEV Protease in 50 mM Tris-HCl pH 7.5, 150 mM NaCl, 1 mM TCEP, 1 mM EDTA, 50 % Glycerol (v/v)

TEV Protease Reaction Buffer

10x conc. reaction buffer containing 500 mM Tris-HCl pH 7.5, 5 mM EDTA, 10 mM DTT

Note:

TEV protease is stable in up to 500 mM NaCl and 300 mM imidazole. Incubation time and temperature may vary depending on the application. TEV protease also cleaves in 16 hours at both 4 °C and 25 °C.