

Recombinant TMPRSS2 Protein, Partial

(Catalog # E80017)

Background

TMPRSS2 plays an important role in the infection mechanism of human coronaviruses, such as SARS-CoV and SARS-CoV-2. Cell entry of human coronaviruses depends on the binding of the viral spike (S) glycoprotein to cellular ACE2 receptor and S protein priming by host cell protease TMPRSS2. This protein is commonly used to study cancer. Previous studies have shown that the encoding gene for TMPRSS2 protein was up-regulated by androgenic hormones in prostate cancer cells and down-regulated in androgen-independent prostate cancer tissue.

Description

The recombinant Transmembrane protease serine 2 (TMPRSS2); expressed in E.coli at the 106-492aa region of TMPRSS2.

Purity

Greater than 90% as determined by SDS-PAGE.

Formulation

Lyophilized powder; Buffer Tris-based buffer, 50% glycerol

Species

Human

Calculated MW

46.9 kDa

Expressed Host

E.coli (Expression Region 106-492aa)

Tag

N-terminal 6xHis-tagged

Uniprot ID

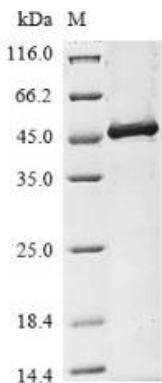
O15393

Storage

Store at -20°C upon receipt, aliquoting is necessary for multiple use. Avoid repeated freeze-thaw cycles.

Alternative Names

D16Ert61e; Epitheliasin; FLJ41954; MGC6821; PP9284; PRSS10; Serine protease 10; TMPRSS2; TMPRSS2 ERG FUSION GENE, INCLUDED; TMPRSS2 ETV1 FUSION GENE, INCLUDED; TMPS2_HUMAN; Transmembrane protease serine 2 catalytic chain; Transmembrane protease, serine 2; Transmembrane protease, serine 2, EC 3.4.219



Tris-Glycine gel: Discontinuous SDS-PAGE (reduced) with 5% enrichment gel and 15% separation gel.