

PRODUCT DATA SHEET



ACE2 (human, recombinant)

CATALOG NO.: SE-520

LOT NO.: XXXXX

DESCRIPTION: MW=85 kDa calculated. Recombinant human angiotensin I converting enzyme-2 (ACE2, ACEH) extracellular domain (residues 1-740) with a C-terminal purification tag, expressed in tissue culture. ACE2 is a cell membrane-inserted zinc carboxypeptidase involved in processing of neuropeptides, possessing 42% sequence identity to ACE1¹⁻³. Drug discovery interests include gastritis and inflammatory bowel disease.

PURITY: >90% by SDS-PAGE. Due to glycosylation, it runs at approximately 100 kDa in SDS-PAGE.

USAGE: Study enzyme kinetics, cleave target substrates, and screen for inhibitors⁴.

SUPPLIED AS: 5µg at 0.25 µg/µl in PBS.

STORAGE: -70°C. The enzyme is stable on ice for at least several hours. However, it is recommended that thawing and dilution of the enzyme be done within as short a time as possible before start of the assay. After initial defrost, aliquot product into individual tubes and refreeze at -70°C. Avoid repeated freeze/defrost cycles.

NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to loss of activity.

REFERENCES:

1. A.J. Turner et al. Can. J. Physiol. Pharmacol. 2002 80 346
2. S.R. Tipnis et al. J. Biol. Chem. 2000 275 33238
3. M. Donoghue et al. Circ. Res. 2000 87 E1
4. L. Huang et al. J. Biol. Chem. 2003 278 15532

FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.