

s-COMT (V108M) (His) (Soluble Catechol-O-Methyltransferase)

CATALOG NO.: MT-11-357

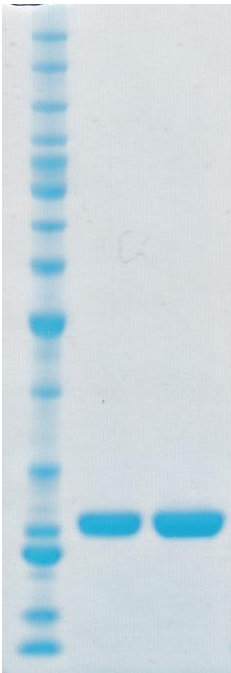
LOT NO.:

DESCRIPTION: Human recombinant s-COMT with methionine (M) in place of valine (V108) (residues 2-221; Genbank Accession # NM_007310; MW = 26.2 kDa) expressed with a C-terminal His tag in *E. coli*.

PURITY: >95% by SDS-PAGE

SUPPLIED AS: µg/µL in 50 mM Tris HCl, pH 7.5, 150 mM NaCl, 1 mM TCEP, 10% glycerol

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 2 and 4 µg of RBC s-COMT (V108M). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, **20**, 15, 10 kDa.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.