

Human

tryptophan

Data Sheet

Description:

<u>TPH2</u>

Human, recombinant, N-terminal His-tag Catalog #: 71193 Lot #: 140902-1 Conc.: 1.08 mg/ml	hydroxylase 2 (TPH2) also known as neuronal tryptophan hydroxylase, GenBank Accession No. NM_173353, a.a.
Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 160 mM	MW = 37 kDa, expressed in an <i>E. coli</i> cell expression system.
imidazole <u>Stability</u> : >6 months at -80 °C. Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 - 0.5%). <u>References</u> : 1. Zill P. et al. Mol. Psychiatr 2004: 9:	Assay Conditions: Reaction was performed at room temperature in assay buffer containing 50 mM MOPS, pH 7.0, 60 μM tryptophan, 100 mM ammonium sulfate, 100 μM ferrous ammonium sulfate, 0.5 mM TCEP, 0.3 mM 6-methyl tetrahydropterin, 0.05 mg/ml catalase, 1 mM DTT and TPH2. Fluorescence signal was measured at Em360nm/Ex300nm.
1030-1036. 2. Sheehan, K., <i>et al., Mol. Psychiatr.</i> 2005; 10 : 944-949.	<u>Application</u> : Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694** Or you can Email us at: <u>info@bpsbioscience.com</u> Please visit our website at: <u>www.bpsbioscience.com</u>