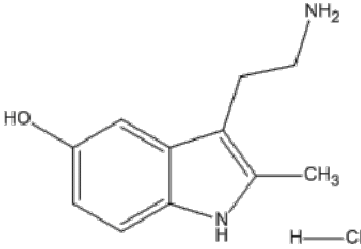


2-Methyl-5-hydroxytryptamine hydrochloride

Instruction Manual

Catalog Number	PK-CA577-2768-25
Description	2-Methyl-5-hydroxytryptamine is a 5-HT ₃ agonist (K _i = 1200 nM) and potent 5-HT ₆ ligand (K _i = 46 nM). Recently, this has been used as one of the components of the chemical cocktails used for reprogramming of mouse somatic cells toward pluripotent cells in the absence of Oct4.
Quantity	25 mg
Sequence / Molecular Weight / Molecular Formula	226.71 Da; C ₁₁ H ₁₄ N ₂ O.HCl
Chemical Structure	
Synonyms	NA
Purity	≥98% as determined by HPLC.
Appearance / Formulation / Solubility	Beige solid. Soluble in water (>20 mg/ml).
Storage & Stability	Store desiccated at RT or 4°C. Protect from light and moisture.
Applications	Stem Cell Fate Regulator. See Description.
References	Hou, P., et al. (2013). Science 341, 651-654.
Caution	Do not take internally. Wear gloves and mask when handling the product! Avoid contact by all modes of exposure.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.