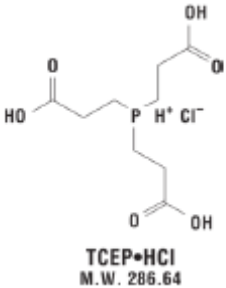


TCEP hydrochloride; (tris(2-carboxyethyl)phosphine)

Instruction Manual

Catalog Number	PK-CA707-91049
Description	TCEP hydrochloride is an odorless reducing agent that is more stable and effective than dithiothreitol (DTT) or 2-Mercaptoethanol. TCEP hydrochloride is soluble in water to 310 grams per liter. Unlike DTT, TCEP hydrochloride retains its reducing power at acid pH (pH 5) and at pH above 7.5. TCEP hydrochloride is unreactive towards other functional groups found in proteins. For most applications, 5-50 mM TCEP can be used to reduce disulfide bonds within a few minutes at room temperature.
Quantity	1 g
Excitation / Emission Maxima	NA
Molecular Structure	 <p style="text-align: center;">TCEP·HCl M.W. 286.64</p>
Molecular Weight / Molecular Formula	286.64 Da; C ₉ H ₁₅ O ₆ P • HCL
Purity	≥98% (as determined by TLC)
Appearance / Formulation / Solubility	White solid. Soluble in water.
Storage & Stability	Store desiccated at -20°C. Protect from light. Keep vial tightly sealed to prevent oxidation.
Applications	see Description
References	1) Anal. Biochem. 180 , 231-236 (1989); 2) Anal Biochem. 220 , 5-10 (1994); 3) Nature Biotech 19 , 379-382 (2001).
Caution	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

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