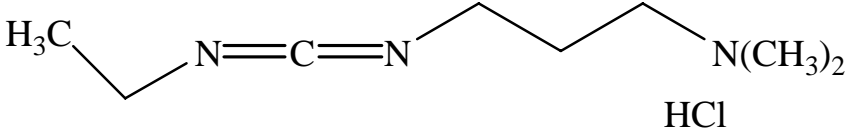


## 1-(3-dimethylaminopropyl-3-ethylcarbodiimide, hydrochloride)

### Instruction Manual

Catalog Number	PK-CA707-59002
Description	EDC (or EDAC) has recently been found to be very useful for fixing chelators <i>in situ</i> , including the fluorescent ion indicators provided by PromoKine. The fixation of ion indicators makes it useful for post histological studies following the physiological experiments. EDC is also a widely used reagent to activate carboxy groups for amine couplings.
Quantity	100 mg
Excitation / Emission Maxima	NA
Molecular Structure	
Molecular Weight / Molecular Formula	191.7 Da; C <sub>8</sub> H <sub>16</sub> ClN <sub>3</sub>
Purity	>98%
Appearance / Formulation / Solubility	White solid; soluble in water, DMF or DMSO.
Storage & Stability	Store desiccated at -20°C.
Applications	See Description
References	1) Tymianski, M. et al. Cell Calcium 21(3), 175(1997) 2) Tetrahedron Lett. 34, 7685(1993)
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.