

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: EDC (EDAC; 1-(3-dimethylaminopropyl-3-ethylcarbodiimide, hydrochloride))

CATALOG #: 59002

MOLECULAR $C_8H_{18}CIN_3$ **INFORMATION:** MW: 191.7

[25952-53-8]

PROPERTIES:

Color & Form White solid **Purity** $\geq 98\%$

Soluble in water, DMSO or DMF

STORAGE AND HANDLING:

Store desiccated at -20 °C.

APPLICATION: EDC (or EDAC) has recently been found to be very useful for fixing chelators *in situ*, including

the fluorescent ion indicators in this catalog. 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.

Ref: Tymianski, M. et al. Cell Calcium 21(3), 175(1997); Tetrahedron Lett. 34, 7685(1993)

TOXICITY: Unknown.

FIRST AID: 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.

Disclaimer: Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.

Revised: April 30, 2018 Page 1 of 1

PSF006