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: CHAPS, ULTRA PURE

CATALOG # 22024-10g, 22024-100g

MOLECULAR $C_{32}H_{58}N_2O_7S$ INFORMATION:MWt: 614.88

[75621-03-3]

PROPERTIES:

Color & FormWhite solidPurity≥98% by HPLCSolubilitySoluble in water

STORAGE AND HANDLING:

Stored desiccated at room temperature.

APPLICATION: CHAPS, 3-[(3-cholamidopropyl)dimethylammonio]-1-propanesulfonate, is a

zwitterionic detergent that is used for protecting the native state of proteins. This zwitterionic detergent is useful for membrane protein solubilization when it is important to maintain protein activity. It is electrically neutral and easily removed by

dialysis.

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.