



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## CHAPS, Ultra Pure

Zwitterionic non-denaturing detergent for solubilizing membrane proteins for protein gels and assays.



### Product attributes

Storage Conditions	Store at room temperature, Desiccate
CAS number	75621-03-3
Molecular weight of antigen	614.88

## Product Description

CHAPS (3-[(3-cholamidopropyl)dimethylammonio]-1-propanesulfonate) is a zwitterionic detergent that is used for protecting the native state of proteins. This detergent is useful for membrane protein solubilization when it is important to maintain protein activity. It is electrically neutral and easily removed by dialysis.

- Zwitterionic detergent
- Solubilizes membranes leaving protein activity intact
- White solid soluble in water
- $C_{32}H_{58}N_2O_7S$

This datasheet was generated on June 12, 2026 at 12:50:57 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/chaps-ultra-pure/>