

Detergent

Catalog No: 37517

Format: 25 ml

Quality Control: Active Motif's Detergent (part number 101683) is a non-ionic, non-denaturing detergent. It can be used for solubilization in protein purification protocols and to increase stringency in wash buffers.

Detergent is provided ready to use.

Storage and Guarantee: Store at room temperature. This product is for research use only and is not for use in diagnostic procedures. This product is guaranteed for 6 months from date of arrival.