

Tagmentase Dilution Buffer

Cat. No. **C01070011**

Format: 200 µl

Product description

Diagenode Tagmentase Dilution Buffer is recommended for dilution of Tagmentase (Cat. No. C01070010) before or after transposome assembly.

Storage conditions

Store at -20°C. Guaranteed stable for 2 years from date of receipt when stored properly.

Storage buffer

Supplied in solution containing 50% v/v glycerol.

Precautions

This product is for research use only. Not for use in diagnostic or therapeutic procedures.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
Rue du Bois Saint-Jean, 3
4102 Seraing - Belgium
Tel: +32 4 364 20 50
Fax: +32 4 364 20 51
orders@diagenode.com
info@diagenode.com

Diagenode LLC. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
Denville, NJ 07834 - USA
Tel: +1 862 209-4680
Fax: +1 862 209-4681
orders.na@diagenode.com
info.na@diagenode.com

Last update: June, 2023