

## Tagmentase Dilution Buffer

**Cat. No.** C01070011

**Lot:** 1/b

**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

This solution contains 50% v/v glycerol.

### Precautions:

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