



## Universal Restriction Buffer

10 x conc.

Cat. No.	Amount
EN-300	5 x 1 ml

### Description:

Universal Restriction Buffer is a single buffer system optimized for fast and efficient digestion fitting to a broad range of restriction enzymes. The proprietary composition ensures highest enzymatic activity without the need for additional optimization steps.

**For general laboratory use.**

**Shipping:** shipped at ambient temperature

**Storage Conditions:** store at 4 °C or -20 °C

**Shelf Life:** 12 months

**Form:** liquid

**Concentration:** 10 x