

CUT&Tag-IT® Assay Buffer Set - Cells

Catalog No: 53169

Format: 1 set

Quality Control: The CUT&Tag-IT™ Assay Buffer Set is useful for performing CUT&Tag reactions using Active Motif's CUT&Tag-IT™ Assembled pA-Tn5 Transposomes (Cat. nos. 53164, 53166).

The CUT&Tag-IT™ Assay Buffer Set includes: Tagmentation Buffer, Dig-Wash Buffer, Dig-300 Buffer, Antibody Buffer, 1X Binding Buffer, and 1X Wash Buffer for use with CUT&Tag-IT™ Assembled pA-Tn5 Transposomes (Cat. Nos. 53164, 53166).

For a complete protocol please refer to the CUT&Tag-IT™ Assay Kit (Cat. Nos. 53160 or 53165) manuals at www.activemotif.com.

Contents:

Tagmentation Buffer, 2 vials of 2 mL, store at -20°C

Dig-Wash Buffer, 50 mL, store at 4°C

Dig-300 Buffer, 50 mL, store at 4°C

Antibody Buffer, 800 µL, store at 4°C

1X Binding Buffer, 55 mL, store at 4°C

1X Wash Buffer, 40 mL, store at 4°C

Storage and Guarantee: Store components according to their labels. This product is guaranteed for 6 months from date of arrival when stored properly.