

ATAC-Seq Buffer Set

Catalog No: 53153

Lot No: 25024183

Format: 1 set

Background: The ATAC-Seq Buffer Set includes ATAC Lysis Buffer, 2X Tagmentation Buffer, and 1% Digitonin for use with Assembled Tn5 Transposomes in 96 ATAC-Seq Assay reactions. We recommend the Pre-indexed Assembled Tn5 Transposomes, 1 x 96 rxns (Cat. No. 53152).

A protocol to use the ATAC-Seq Buffer Set in ATAC-Seq is included on the Assembled Tn5 Transposomes (Cat. No. 53152) web page.

Contents:

ATAC Lysis Buffer, 15 mL, store at RT or 4°C

2X Tagmentation Buffer, 2 vials of 1.5 mL, store at -20°C

1% Digitonin, 2 vials of 25 µL, store at -20°C

Quality Control: The ATAC-Seq Buffer Set is useful for performing ATAC-Seq reactions using Active Motif's Pre-indexed and Assembled Tn5 Transposomes (Cat. No. 53152) or with Tn5 (Cat. Nos. 81284 or 81286) that has been "loaded" or assembled into transposomes.

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