

Q5 Polymerase and Library Purification Kit

Catalog No: 53159

Format: 1 set

Background: The Q5 Polymerase and Library Purification Kit includes:

5X Q5 Reaction Buffer, 960 μ L, store at -20°C ,
dNTP Mix, 10 mM, 96 μ L, store at -20°C ,
Q5 DNA Polymerase, 48 μ L store at -20°C ,
SPRI Beads, 5.76 mL store at 4°C ,
Silica Beads, 288 μ L store at 4°C ,
DNA Purification Binding Buffer, 4.8 mL store at RT,
DNA Purification Elution Buffer, 5.28 mL store at RT,
DNA Purification Wash Buffer, 7.68 mL store at RT.

Quality Control: The Q5 Polymerase and Library Purification Kit is useful for DNA amplification and library purification in ATAC-Seq reactions in High-Throughput ATAC-Seq Express Kit (Cat. No. 53158) or other Tn5-based applications.

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