

PRODUCT NAME

Blocker for ChIP-Ab binding beads

Cat. No: kch-505-200

Format: 200 µl

Cat. No: kch-505-500

Format: 500 µl

Description

This reagent is a BSA solution (10x) which has been extensively validated in chromatin immunoprecipitation (ChIP) assays. It is intended for being used as a blocker for ChIP-Ab binding beads (agarose and magnetic).

Format

Supplied as a 5% solution.

Storage

Store at -20°C. Stable for 2 years.

Precautions

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