

CUTANA™ Bead Activation Buffer

Catalog No	21-1001	Pack Size	48 Reactions
Lot No	24165001-81	Applications	CUT&RUN, CUT&Tag

DESCRIPTION

CUTANA™ Bead Activation Buffer is used to prepare Concanavalin A (ConA) conjugated paramagnetic beads for CUT&RUN and CUT&Tag. Activated ConA beads are used in both CUT&RUN and CUT&Tag workflows to immobilize cells or nuclei.

TECHNICAL INFORMATION

Storage	Stable for 12 months at 4°C from date of receipt.
Formulation	20 mM HEPES pH 7.9, 10 mM KCl, 1 mM CaCl ₂ , 1 mM MnCl ₂ .
Instructions for Use	CUTANA™ Bead Activation Buffer includes sufficient buffer for 48 CUT&RUN or CUT&Tag reactions using the CUTANA™ protocols (www.epicypher.com/protocols).

RECOMMENDED COMPANION PRODUCTS

Item	Catalog No.
CUTANA™ CUT&RUN Kit	14-1048 / 14-1048-24
CUTANA™ CUT&RUN Library Prep Kit	14-1001 / 14-1002
CUTANA™ CUT&Tag Kit	14-1102 / 14-1103
CUTANA™ pAG-MNase	15-1016 / 15-1116
CUTANA™ pAG-Tn5	15-1017 / 15-1117
CUTANA™ ConA Beads	21-1401 / 21-1411
CUTANA™ Nuclei Extraction Buffer	21-1026
CUTANA™ Stop Buffer	21-1003
CUTANA™ 5% Digitonin	21-1004
CUTANA™ E. coli Spike-in DNA	18-1401
Magnetic Separation Rack (0.2 mL / 1.5 mL tubes)	10-0008 / 10-0012
8-strip 0.2 mL Tubes	10-0009

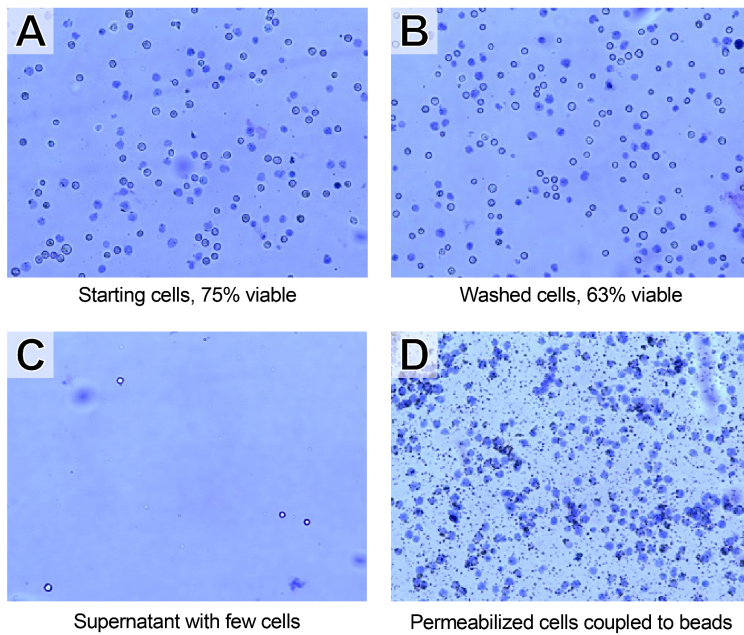


FIGURE 1 Cell binding to activated ConA beads. ConA beads were washed with CUTANA™ Bead Activation Buffer and loaded with K562 cells as described in the CUTANA™ CUT&RUN and CUT&Tag protocols (www.epicypher.com/protocols). Trypan Blue staining shows cells before (panels A and B) and after (C and D) bead binding. After incubation with activated beads, the supernatant contains few unbound cells (C), demonstrating efficient binding. Permeabilized cells bound to activated beads are observed in the cell-bead slurry (D).