

NCOA3/SRC3 CUTANA™ CUT&RUN Antibody



EpiCypher®

Catalog No. 13-2013

Lot No. 21085001-04

Pack Size 100 µL

Type Polyclonal **Target Size** 155 kDa

Host Rabbit **Format** Aff. Pur. IgG

Reactivity Human

Applications CUT&RUN, WB

Product Description:

This antibody meets EpiCypher's "CUTANA Compatible" criteria for performance in Cleavage Under Targets and Release Using Nuclease (CUT&RUN) and/or Cleavage Under Targets and Tagmentation (CUT&Tag) approaches to genomic mapping. Every lot of a CUTANA Compatible antibody is tested in the indicated CUTANA approach using EpiCypher optimized protocols (EpiCypher.com/resources/protocols) and determined to yield peaks that show a genomic distribution pattern consistent with reported function(s) of the target protein. Consistent with its role as a nuclear receptor coactivator that facilitates ER-mediated transcription (1), NCOA3 (SRC3) antibody shows CUT&RUN peaks in response to estradiol stimulation (**Figure 1**) that overlap with known estrogen response element (ERE) binding motifs (**Figure 2**). SRC3 CUT&RUN peaks also overlap with ER alpha antibodies (N- and C-term) (**Figure 2**).

Immunogen:

A synthetic peptide corresponding to human NCOA3/SRC3 amino acids 900 - 950.

Formulation:

Antigen affinity-purified antibody (1.0 mg/mL) in Tris-citrate/phosphate buffer pH 7 to 8, 0.09% sodium azide.

Storage and Stability:

Stable for 1 year at 4°C from date of receipt.

Application Notes:

Recommended Dilutions:

CUT&RUN: 0.5 µg

WB: 1:2,000 - 1:10,000

References:

1. Wagner et al. (2013) *BMC Cancer* 4:13-570

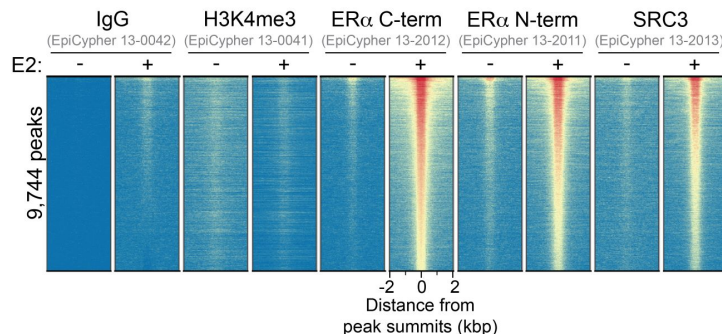


Figure 1: SRC3 enrichment in CUT&RUN. Serum-starved MCF7 cells were treated with 100 nM estradiol (E2) or vehicle control for 45 minutes. CUT&RUN was performed using 500,000 cells with 0.5 µg of each indicated antibody (gray text). Heatmap shows CUT&RUN enrichment in aligned rows ranked by intensity (top to bottom; relative to ER alpha C-term). Red indicates high localized enrichment and blue denotes background signal.

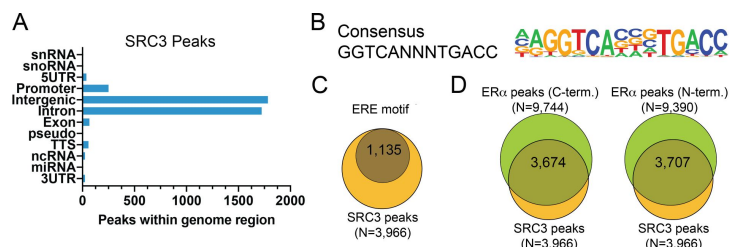


Figure 2: SRC3 peak analysis in CUT&RUN. Peaks from the E2-treated samples in Figure 1 were called using MACS2. (A) The number of SRC3 peaks which fall into distinct classes of functionally annotated genomic regions is plotted. (B) Homer analysis determined that the ERE consensus motif, represented as a sequence logo position weight matrix, was enriched under SRC3 peaks. (C) The number of SRC3 peaks containing consensus motifs from panel B is shown by Venn Diagram. (D) The number of SRC3 peaks that overlap with ER alpha C-term and ER alpha N-term antibodies are represented by Venn Diagram.

This product is provided for commercial sale under license from Bethyl Laboratories, Inc.

Applications Key: ChIP: Chromatin immunoprecipitation; CUT&RUN: Cleavage Under Targets and Release Using Nuclease; CUT&Tag: Cleavage Under Targets and Tagmentation; E: ELISA; FACS: Flow cytometry; ICC: Immunocytochemistry; IF: Immunofluorescence; IHC: Immunohistochemistry; IP: Immunoprecipitation; L: Luminex; WB: Western Blot. **Reactivity Key:** B: Bovine; Ce: *C. elegans*; Ch: Chicken; Dm: *Drosophila*; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: *S. cerevisiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish

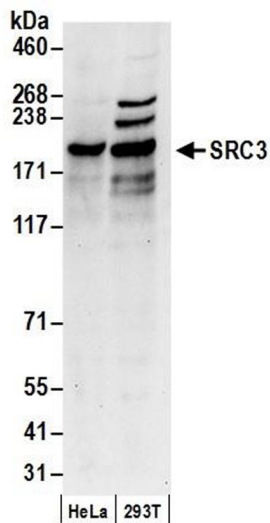


Figure 3: Western blot detection of human SRC3. Whole cell lysates were isolated from HeLa and HEK293T cells using NETN lysis buffer. Lysate (50 μ g) was loaded onto 4-8% SDS-PAGE gel and analyzed under standard western blot conditions using SRC3 antibody (0.1 μ g/mL).

This product is provided for commercial sale under license from Bethyl Laboratories, Inc.

Applications Key: ChIP: Chromatin immunoprecipitation; CUT&RUN: Cleavage Under Targets and Release Using Nuclease; CUT&Tag: Cleavage Under Targets and Tagmentation; E: ELISA; FACS: Flow cytometry; ICC: Immunocytochemistry; IF: Immunofluorescence; IHC: Immunohistochemistry; IP: Immunoprecipitation; L: Luminex; WB: Western Blot. **Reactivity Key:** B: Bovine; Ce: *C. elegans*; Ch: Chicken; Dm: *Drosophila*; Eu: Eukaryote; H: Human; M: Mouse; Ma: Mammal; R: Rat; Sc: *S.cerevesiae*; Sp: *S. pombe*; WR: Wide Range (predicted); X: Xenopus; Z: Zebrafish