

ChIL Probe (goat, anti-rabbit)

Cat. No. C01060020

Format: 12 µl (24 rxns)

Suggested dilutions: 1:100-1:200 in 0.5M NaCl/buffer of your choice

Product specification:

Antibody: AffiniPure Goat Anti-Rabbit IgG (H+L)

Target: Rabbit

Host: Goat

Antibody Format: Whole IgG

Specificity: IgG (H+L)

Minimal Cross Reactivity: Human, Mouse, Rat Serum Proteins

Conjugate: Unconjugated

Product Category: Whole IgG Affinity-Purified Antibodies

Clonality: Polyclonal

Conjugation Compound: DBCO-PEG5-NHS Ester, MW = 693.74 [Click Chemistry Tools cat#: A102P-2]

ChIL oligo DNA:

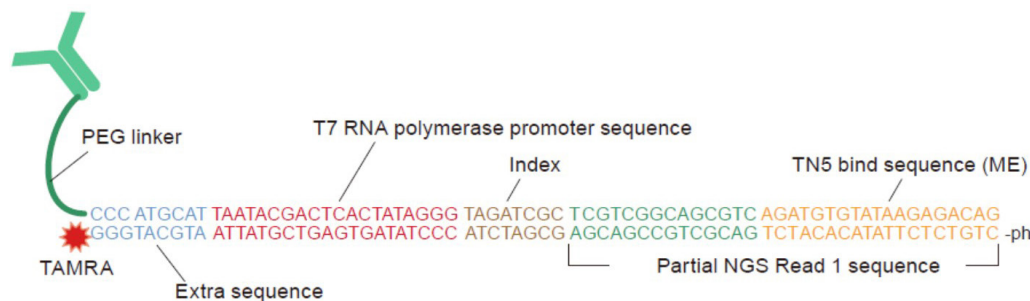


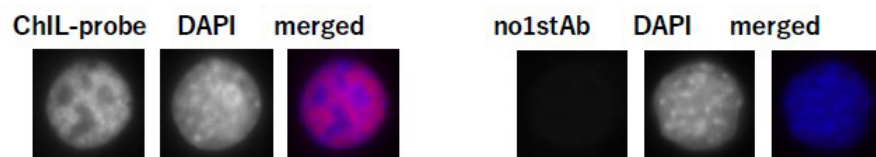
Figure 1. Structure of ChIL probe

Storage conditions: Store at -20°C under sterile/dark condition. Prepare working dilution on day of use.

Storage buffer: Supplied in Phosphate Buffered Saline (pH 7.4)

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

VALIDATION DATA



(exposure time : 1/25 sec all)

Figure 2. Staining pattern of ChIL probe

Primary antibody: H3K4me3

Cell: C2C12

ChIL probe Staining buffer: 0.5M NaCl/0.1xblocking one P(#05999-84)/PBS

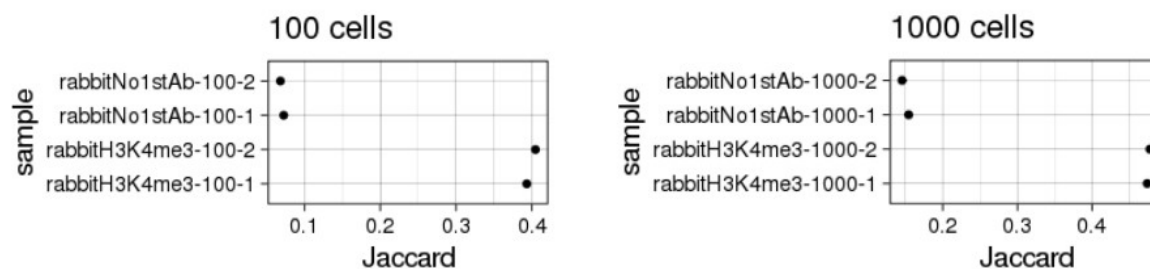


Figure 3. Plotting Jaccard similarity index

ChIL data(H3K4me3) vs ENCODE ChIP-Seq (H3K4me3). Higher number indicates two data sets are more similar.

No1stAb: no primary antibody + ChIL probe. (= negative control)

Reference:

Chromatin integration labeling for mapping DNA-binding proteins and modifications with low input.

Handa T#, Harada A#, Maehara K#, Sato S, Nakao M, Goto N, Kurumizaka H, Ohkawa Y*, Kimura H**Nat Protoc.* 2020