

ETO polyclonal antibody

Other names: RUNX1T1, AML1T1, CBFA2T1, CDR, MTG8, ZMYND2

Cat. No. C15310001

Type: Polyclonal

ChIP-grade / ChIP-seq-grade

Source: Rabbit

Lot #: A710-001

Size: 100 µl

Concentration: not determined

Specificity: Human: positive

Other species: not tested

Purity: Whole antiserum from rabbit containing 0.05% azide.

Storage: Store at -20°C; for long storage, store at -80°C.

Avoid multiple freeze-thaw cycles.

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

Description: Polyclonal antibody raised in rabbit against human ETO (runt-related transcription factor 1; translocated to, 1 (cyclin D-related)) using two KLH-conjugated synthetic peptides containing sequences from the N-terminal and the central region of the protein, respectively.

Applications

	Suggested dilution	Results
ChIP*	4 µl/ChIP	Fig 1, 2
ELISA	1:100	Fig 3

* Please note that the optimal antibody amount per ChIP should be determined by the end-user. We recommend testing 1-10 µl per IP.

References citing this antibody:

(1) Martens JHA, Mandoli A, Simmer F, Wierenga B-J, Saeed S, Singh AA, Altucci L, Vellenga E, Stunnenberg HG (2012) ERG and FLI1 binding sites demarcate targets for aberrant epigenetic regulation by AML1-ETO in acute myeloid leukemia. Blood 120: 4038-4048.

Target description

ETO (UniProtKB/Swiss-Prot entry Q06455) is a transcriptional regulator which belongs to the myeloid translocation gene family. ETO exerts its function by interaction with transcription factors bound to promoters and binding to histone deacetylases. It recruits a range of corepressors to facilitate transcriptional repression. The t(8;21)(q22;q22) translocation is one of the most frequent karyotypic abnormalities in acute myeloid leukaemia. This translocation produces a chimeric gene made up of the 5'-region of AML1 and the 3'-region of the ETO gene. The chimeric protein is thought to associate with the nuclear corepressor/histone deacetylase complex to block hematopoietic differentiation.

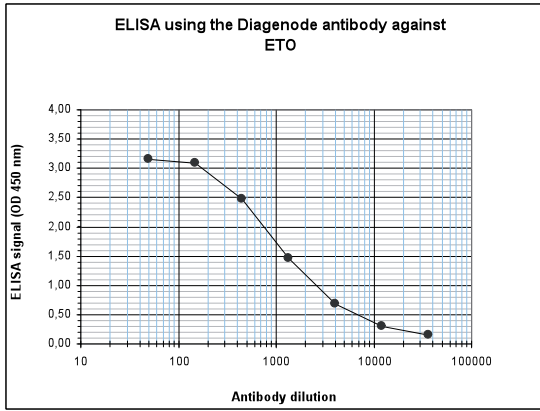


Figure 3. Determination of the antibody titer

To determine the titer of the antibody, an ELISA was performed using a serial dilution of the Diagenode antibody directed against human ETO (Cat. No. C15310001). The plates were coated with the peptides used for immunization of the rabbit. By plotting the absorbance against the antibody dilution (Figure 3), the titer of the antibody was estimated to be 1:1,300.

Diagenode sa. BELGIUM | EUROPE

LIEGE SCIENCE PARK
 Rue Bois Saint-Jean, 3
 4102 Seraing (Ougrée) - Belgium
 Tel: +32 4 364 20 50
 Fax: +32 4 364 20 51
 orders@diagenode.com
 info@diagenode.com

Diagenode Inc. USA | NORTH AMERICA

400 Morris Avenue, Suite 101
 Denville, NJ 07834 - USA
 Tel: +1 862 209-4680
 Fax: +1 862 209-4681
 orders.na@diagenode.com
 info.na@diagenode.com

Last update: August 31, 2012