

Antibody Package (anti-rabbit)

Cat. No. C01070022, C01070026

Format: 24 rxns, 8 rxns

Product content

Antibody Package	Volume – 8 rxns Cat. No. C01070026	Volume – 24 rxns Cat. No. C01070022	Storage
ChIP-seq grade rabbit antibody H3K27me3 (1 µg/µl)	2 µl	2 µl	-20°C
Rabbit IgG (1 µg/µl)	2 µl	2 µl	-20°C
Anti-rabbit Secondary antibody	8 µl	24 µl	-20°C
ChIP-seq grade TSH2B primer pair (human)*	8 µl	24 µl	-20°C
ChIP-seq grade GAPDH TSS primer pair (human)*	8 µl	24 µl	-20°C

Product description

The Antibody Package (anti-rabbit), Cat. No. C01070022, is designed for use across chromatin profiling applications and is compatible with assays such as ChIP and CUT&Tag. This package is recommended when the primary antibody of interest used in the assay has been raised in rabbit.

The Antibody Package (anti-rabbit) contains species-specific controls antibodies (anti-rabbit secondary antibody, rabbit H3K27me3 positive control antibody, rabbit negative IgG control) and primer pairs (TSH2B and GAPDH TSS) for the evaluation of relative enrichment by qPCR. The TSH2B and GAPDH TSS primer pairs are positive and negative primers, respectively, for H3K27me3 control antibody and are only compatible with human samples.

The package contains enough secondary antibody to perform 24 rxns. The positive and negative control antibodies are enough for 2 reactions each. ChIP-seq grade TSH2B (positive) and GAPDH TSS (negative) primer pairs are enough for 24 reactions each.

Diagenode, SA
BELGIUM | EUROPE

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

Diagenode, LLC
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: June, 2026