

EZH2 phospho Thr345 antibody (pAb)

Catalog Nos: 61241, 61242**RRID:** AB_2793564**Isotype:** IgG**Application(s):** WB**Reactivity:** Human, Mouse**Volumes:** 100 µl, 10 µl**Purification:** Affinity Purified**Host:** Rabbit**Molecular Weight:** 96 kDa

Background: **EZH2** – (Enhancer of Zeste homolog 2, also designated Enx1) is a human homolog of the Drosophila Polycomb-group protein Enhancer of Zeste protein. It contains a SET domain that catalyzes the methylation of histone H3 at lysine 27 (Histone H3 Lys27 monomethylated, dimethylated and trimethylated). Polycomb-group proteins repress gene expression by binding to chromatin and locally altering chromatin structure. **EZH2**, BMI-1 and Suz12 are present in the PRC2 and PRC3 protein complexes that function as mediators of epigenetic transcriptional silencing. Threonine 345 of mouse Ezh2 is phosphorylated in a cell cycle-dependent manner and may play a role in the interaction of Ezh2 with HOTAIR and the 59-RepA repeats of Xist. (see reference)

Immunogen: This EZH2 phospho Thr345 antibody was raised against a peptide containing phosphothreonine 345 of mouse EZH2 (enhancer of zeste homolog 2 (Drosophila))

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

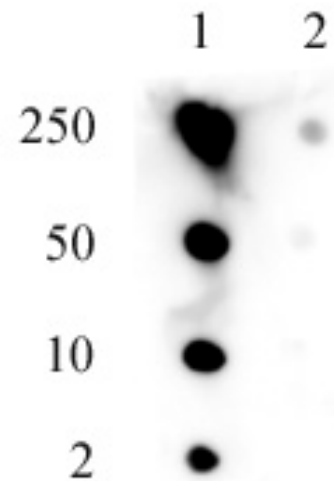
Published Applications:

WB

See references for more information. Individual optimization may be required.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.

**EZH2 phospho Thr345 pAb tested by dot blot analysis.**

Dot blot analysis was used to confirm the specificity of EZH2 phospho Thr345 antibody. Modified and unmodified peptide were spotted onto PVDF and probed with the antibody at a dilution of 1:5,000. Decreasing amounts of peptide were spotted in each row. Lane 1: Peptide phosphorylated at threonine 345. Lane 2: Unmodified EZH2 peptide.