

Anti-conjugated L-Serine antibody

Ref: IS1003

| | |
|-------------------|---------------------------------|
| Clonality | Polyclonal |
| Host | Rabbit |
| Applications | <u>IHC</u> / IF |
| Specie reactivity | Reacts with all species |
| References | Not yet cited to our knowledge. |
| Format | 50µl |

Product information

Product overview

| | |
|---------------------|--|
| Product name | Anti-conjugated L-Serine antibody |
| Synonyms | Anti-L-Serine acid antibody Anti-(S)-2-amino-3-hydroxypropionic acid antibody |
| Immunogen | Conjugated L-Serine |
| Specificity | When tested in competitive ELISA, the anti-conjugated L-Serine antibody did not show any significant cross reactivity with L-Serine analogs, including D-Serine and L-Threonine conjugates |
| Lot number | 140301 |

Reconstitution & storage

| | |
|-----------------------|--|
| Form | Lyophilized powder |
| Purity | Purified anti-serum |
| Storage | Store at 4 °C |
| Storage buffer | Before use, vial should be resuspended in 50 µL of ultrapure water. Store at +4 °C for short term (1-2 weeks). Aliquot and store at -20 °C for long term. Avoid repeated freeze / thaw cycles. |

Applications

| | |
|-----------------------------------|--|
| Immunohistochemistry (IHC) | Optimal working conditions must be determined by the end-user. |
| Immunofluorescence (IF) | Optimal working conditions must be determined by the end-user. |
| Restrictions | For research use only |

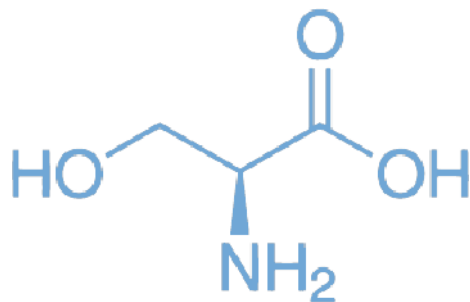
References

Selected publications about L-Serine:

- [Pollari S et al. Enhanced serine production by bone metastatic breast cancer cells stimulates osteoclastogenesis. Breast Cancer Res Treat. 2011 Jan;125\(2\):421-30. doi: 10.1007/s10549-010-0848-5. Epub 2010 Mar 30.](#)
- [Yang JH et al. Brain-specific Phgdh deletion reveals a pivotal role for L-serine biosynthesis in controlling the level of D-serine, an N-methyl-D-aspartate receptor co-agonist, in adult brain. J Biol Chem. 2010 Dec 31;285\(53\):41380-90. doi: 10.1074/jbc.M110.187443. Epub 2010 Oct 21.](#)
- [Tabatabaie L, Klomp LW, Berger R, de Koning TJ. L-serine synthesis in the central nervous system: a review on serine deficiency disorders. Mol Genet Metab. 2010 Mar;99\(3\):256-62. doi: 10.1016/j.ymgme.2009.10.012. Epub 2009 Oct 25.](#)

Pictures

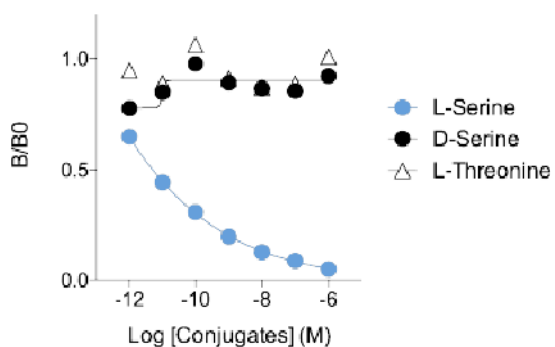
L-Serine



L-Serine

Affinity & specificity of anti-L-Serine antibody

Anti-L-Serine pAb (Rabbit)



Competitive ELISA demonstrates that low amounts of L-Serine conjugate are required to abolish antigen-antibody reaction (high affinity), while rising concentrations of D-Serine and L-Threonine conjugates do not affect reaction (high specificity).

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com • info@eaglebio.com



EAGLE
BIOSCIENCES

For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.