

FHM vitellogenin standard

Contents

Each vial normally contains 3-5 µg purified Fathead Minnow vitellogenin (Vtg). The content of the vial will vary from batch to batch, see exact amount on the label of each vial.

Source of vitellogenin

Fathead Minnow (*Pimephales promelas*) induced with 17β-estradiol.

Purification procedure

FHM Vtg was purified from plasma of 17β-estradiol-induced fish by ion exchange chromatography and gel filtration (1).

Applications

The lyophilized Fathead Minnow Vtg may be used as a positive control in western blot and ELISA. Freshly reconstituted Vtg may also be used as standard in a quantitative ELISA.

Storage

Lyophilized vitellogenin can be stored at 4°C. We recommend reconstitution in 300-1000 µl cold PBS immediately before use. Do not freeze and thaw if Vtg is used as a quantitative standard. For use only as a positive control the solution may be aliquoted and stored at -20°C. Avoid repeated freezing and thawing.

Note: If the solution of vitellogenin after reconstitution appears turbid, add 1-2 µl of 0.2 M EDTA, pH 7.7 until the solution becomes clear.

For research use only

References

1) Brion, F., Nilsen, B.M, Eidem, J.K., Goksøyr, A. and Porcher, J-M. (2002) *Environ Toxicol Chem.* 28, 1699-1708