

Rainbow trout vitellogenin standard

Contents

Each vial normally contains 5-10 µg purified rainbow trout 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

Rainbow trout (*Oncorhynchus mykiss*) induced with 17β-estradiol.

Purification procedure

Rainbow trout Vtg was purified from plasma of 17β-estradiol-induced fish by selective precipitation with MgCl₂ in the presence of EDTA, essentially as described by Norberg and Haux (1) and Arukwe et al (2).

Applications

The lyophilized rainbow trout 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 Vtg 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 Vtg after reconstitution appears turbid, add 1-2 µl of 0.2 M EDTA, pH 7.7 until the solution becomes clear.

References

- 1) Norberg, B. and Haux, C. (1988) Fish Physiol. Biochem. 5, 59-68.
- 2) Arukwe, A., Knudsen, F. R. and Goksøyr, A. (1997) Environ. Health Perspect. 105, 418-422

For research use only