

## Medaka vitellogenin standard

### Contents

Each vial normally contains 5-10 µg purified medaka 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

Japanese medaka (*Oryzias latipes*) induced with 17β-estradiol.

### Purification procedure

Medaka Vtg was purified from whole body homogenate of 17β-estradiol-induced fish by ion exchange chromatography and gel filtration (1).

### Applications

The lyophilized medaka 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.

### References

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

For research use only