

IMCO Corporation Ltd AB

PO Box 21195

SE -100 31 Stockholm

Sweden

www.imcocorp.se

PRODUCT SPECIFICATION

Product No.	ATRX-01
Product description:	Sheep Antiserum to <i>E. coli</i> Thioredoxin 1-(Trx1)
Lot No.	04
Specifications:	Gamma globulin fraction precipitated with (NH ₄) ₂ SO ₄ , dialysed against 0.15 M NaCl, 50 mM Tris-HCl, 1mM EDTA, pH 7.5 and sterile filtered. Volume: 1ml. Sodium azide has been added to a final concentration of 15 mM (0.1%).
Reference:	Randahl, H. et al. (1982) Eur J Biochem, 128, 445-449.
Storage:	Keep at +4°C
Examples of use:	1 ml of this solution will completely inhibit the activity of 12 µM thioredoxin in an assay with thioredoxin reductase, see Holmgren A. and Björnstedt, M. (1995) Methods Enzymol, 252, 199-202.

NOTE! During shipment, the contents of the antibody ampoule may have been redistributed on the inner walls. Before opening the screw-cap of the antibody tube, we recommend to centrifuge the tube shortly, in order to recover liquid from tube cap and walls.