PROTEOSTAT[®] Thioredoxin-1 assay kit

The PROTEOSTAT® Thioredoxin-1 assay kit provides a simple, homogenous assay for screening modulators of Trx-1 enzymatic activity in microplates. This is accomplished by monitoring the Trx-1 catalyzed reduction of insulin in the presence of Dithiothreitol (DTT), resulting in the formation of insulin aggregates which then bind avidly to the red-emitting fluorogenic PROTEOSTAT® Trx-1 detection dye. The fluorescence-based assay provides a vastly improved assay signal window, improved lower detection limit, and superior Z'-score (>0.8) relative to turbidity measurements. Intra-plate and inter-plate CVs using this assay are typically 3-6%. Human recombinant thioredoxin-1 enzyme and the thioredoxin inhibitor, aurothiomalate, are included in the kit, as well as all reagents necessary for monitoring changes in Trx-1 activity. The PROTEOSTAT® Thioredoxin-1 assay kit is capable of providing a quantitative readout of Trx-1 enzymatic activity in a robust and highthroughput fashion and can be applied to identification of Trx-1 inhibitors from chemical libraries.

Citations: 5

View Online »

Ordering Information

Order Online »

ENZ-51033-KP002

2x96 wells

Manuals, SDS & CofA

View Online »

Handling & Storage

Use/Stability Upon receipt, store the Thioredoxin-1 at -80°C. Store the remaing reagents at ≤-20°C

protected from light.

Handling Protect from light. Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Fluorescent detection

Application NotesThis kit provides a simple, homogenous assay for

screening modulators of Thioredoxin-1 enzymatic activity.

Contents PROTEOSTAT[®] PDI detection reagent: 20 μL

Thioredoxin-1 (human) (recombinant): 2 vials x 80 µL

Insulin: 2 vials x 1.8 µmol

Aurothiomalate (inhibitor): 2 vials x 25 nmol

PBE buffer: 25 mL Stop reagent: 1 mL DTT: 2 vials x 1.3 mL Deionized water: 5 mL

Quality Control Using the procedure described in the manual, a sample

from each lot of PROTEOSTAT[®] Thioredoxin-1 assay kit is used to assay the following: (1) Insulin with Trx-1; (2) Insulin without Trx-1; and (3) Insulin with Trx-1 and Aurothiomalate. The Z'-factor is greater than 0.8. Trx-1 is

inhibited by at least 50% with Aurothiomalate.

UniProt ID P10599

Last modified: May 29, 2024



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com