

PRODUCT INFORMATION

DNA/RNA Oxidative Damage Express ELISA AChE Tracer

Item No. 401156



Overview

Contents: This vial contains lyophilized DNA/RNA Oxidative Damage Express ELISA AChE Tracer.
Storage: -20°C
Stability: ≥2 years

Laboratory Procedures

DNA/RNA Oxidative Damage Express ELISA AChE Tracer has been tested and formulated to work exclusively with Cayman's DNA/RNA Oxidative Damage Express ELISA Kit (Item No. 502890). Please visit www.caymanchem.com/catalog/502890 for the kit protocol, procedures, and product handling.

Precaution

This product may not perform as described if used with other assay reagents or protocols.

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/02/2026

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM