

LANCE® *Ultra*

Eu-W1024 labeled Anti-GST Antibody

Product number:	AD0253	Lot Number:	3259627
------------------------	--------	--------------------	---------

Product Format:	AD0252: 10 µg
	AD0253: 50 µg
	AD0254: 1 mg

Manufacturing date:	2/13/2024	Document version:	1
----------------------------	-----------	--------------------------	---

Product Information

Antibody:	The GST antibody is a goat polyclonal antibody that recognizes Glutathione S-Transferase [Schistosoma japonicum].
Storage Buffer:	50 mM Tris-HCl pH 7.4, 150 mM NaCl, 0.05% NaN ₃ , 0.1% BSA
Molecular Weight:	160 000
Stability:	This product is stable for at least 24 months from the manufacturing date when stored in its original packaging and the recommended storage conditions.
Storage Conditions:	Store at 4°C

Quality Control

The QC release specifications are based on spectrophotometric analysis of the labeled antibody. We certify that results meet our quality release criteria.

Labeling Ratio:	6.3 / 1
Concentration:	0.625 µM (100 µg/mL)

Please visit our website for additional information on LANCE *Ultra* technology at www.revvity.com

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

AD0272-3-4-R Rev01