

PhenoVue™ Research Reagents

PhenoVue Fluor 647 - Goat Anti-Rat Antibody Highly Cross-Adsorbed

Product number: 2GXRT647H1

Lot Number: 01A

Appearance: Blue lyophilizate

Quantity: 1 mg

Buffer : 100 mM PO4 + 1 % BSA, pH 7

Expiration date: 13-Sept-2026

Document version: 1

This product has met the quality control standards defined by Revvity, Inc. for cellular imaging application.

	Criteria	Lot data	Specifications
PHOTOPHYSICS	Absorption maximum	Meets specifications	650 ± 5 nm
MICROSCOPY ¹	Background Staining	Meets specifications. Meets specifications	Negligible Good
CROSS-REACTIVITY ²	Species IgG Isotype IgG	Meets specifications. Meets specifications	Conforms to Product Information sheet. Recognises all isotypes

1. Performed with an internal protocol. New batch is compared with the previous one to ensure batch to batch reproducibility.
2. Performed with a FLISA internal protocol on rat, mouse, goat, rabbit, guinea pig, chicken, human, bovine, donkey, hamster, sheep, and horse species IgG and on Rat IgG1, IgG2a, IgG2b, IgG2c Isotypes.

Quality control date	Supervisor's initials
22-09-2023	TR

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.