

Lot specific information

PhenoVue™ Fluor 568 goat anti-Rabbit antibody Highly Cross-adsorbed

Product number: 2GXRB568H1 Lot Number: 01A

Expiration date: Extended to Document version: 1
02/12/2025

Appearance : Purple Lyophilizate
Quantity : 1mg
Buffer : 100mM PO4 + 1% BSA, PH 7

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

	LOT DATA	SPECIFICATION
PHOTOPHYSICS		
Absorption maximum	Meets specifications	578 ± 5nm
MICROSCOPY ¹		
Background	Meets specifications	Negligible
Staining	Meets specifications	Good
CROSS-REACTIVITY ²		
Species IgG	Meets specifications	Conforms to Product Information Sheet
Isotypes IgG	Meets specifications	Recognizes 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 Mouse, Rat, Goat, Rabbit, Guinea Pig, Chicken, Human, Bovine, Donkey, Hamster, Sheep and Horse species IgG and on Mouse IgG1, IgG2a, IgG2b, IgG2c, IgG3 Isotypes.

Quality control date	Supervisor's initials
25/07/2023	LN

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.