

Lot specific information

PhenoVue™ Fluor 488 goat anti-Rabbit antibody Cross-adsorbed

Product number: 2GXRB488C1 Lot Number: 01B

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

The expiration date of batch 01B has been extended to 2025-12-02

Appearance : Orange 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 | 495 ± 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 |
|----------------------|-----------------------|
| 2025/03/20 | ML |

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.

