

**B**

LANCE® Ultra

## Anti-FLAG ULight

<b>Product number:</b>	TRF0059D	<b>Lot Number:</b>	3324078
------------------------	----------	--------------------	---------

---

<b>Product Format:</b>	TRF0059D: 1 nmoles (2 x 1mL)
	TRF0059M 10 nmoles (2 x10mL)
	TRF0059R: 100 nmoles (2 x 100mL)

<b>Manufacturing date:</b>	9/03/2024	<b>Document version:</b>	1
----------------------------	-----------	--------------------------	---

### Product Information

**Application:** ULight has been conjugated to anti-FLAG antibody. The mouse monoclonal recognizes the FLAG sequence at the N-terminus, Met N-terminus, C-terminus, and internal sites. This toolbox can be used to capture FLAG-tagged proteins and peptides,

**Stability:** This product is stable for at least 12 months from the manufacturing date when stored in its original packaging and at the recommended storage conditions.

**Storage Conditions:** Store at 4°C in the dark

### Quality Control

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

<b>Labeling Ratio:</b>	4.9 / 1
<b>Concentration:</b>	0.5 µM , 80 µg/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.

TRF0059-R Rev01