

AlphaLISA®

Lysis Buffer (5X)

Product number:	AL003 C/F	Lot Number:	3316049 , 3296411
------------------------	-----------	--------------------	-------------------

Product Format:	AL003C: 10 mL
	AL003F: 100 mL

Manufacturing date:	8/02/2024	Document version:	1
----------------------------	-----------	--------------------------	---

Product Information

Description: AlphaLISA Lysis Buffer (5X)

Presentation: Liquid, 5X

Storage: Store the 5X solution at 4°C.

Stability: This product is stable for at least 12 months from the manufacturing date when stored in its original packaging at the recommended storage temperature. 1X Buffer supplemented with Protease inhibitor cocktail is stable for 1 day at 4°C. Long-term storage of 1X solution is not recommended. The color of the solution may change over time but is not expected to impact buffer performance.

Reconstitution: Add 4 volumes of HPLC grade H₂O to 1 volume of 5X AlphaLISA Lysis Buffer. Standard protease inhibitor cocktail is recommended to be added prior to use: add 0.25 mL of Protease inhibitor cocktail (Sigma-Aldrich, Cat. No. P2714, or equivalent; reconstituted according to the manufacturer's instructions) per 5 mL of 1X lysis buffer.

Please visit our website for additional information on AlphaLISA 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.

AL003-R Rev 01