

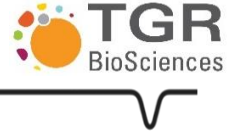
AlphaLISA® SureFire® Ultra™

AlphaLISA CaptSure Acceptor Beads (2mg/ml)

Part numbers: ALSU-ACAB-1.2ML

Lot Number: U17836

Manufacturing date: July 16, 2024

**CONTENTS:****Batch #:**

AlphaLISA CaptSure Acceptor Beads (2mg/ml) Q2147-CE

STORAGE / HANDLING:

Store all Buffers at 4°C

Donor beads are used under low light conditions and should be stored in the box at 4°C. Vortex Acceptor and Donor Beads vigorously prior to taking samples for use.

KIT STABILITY:Stable for at least **12** months from the manufacturing date when stored under recommended storage conditions.**This product is not for resale or distribution except by authorized distributors.**For more information on protocols, please refer to the generic AlphaLISA *SureFire Ultra* Assay Kits Manual, available from 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.