

Caution: For Laboratory Use. A research chemical for research purposes only

AlphaLISA MES Buffer (5X)

Product No.: AL017C/F

Product Information

- Application:** The AlphaLISA MES buffer (5X) has been especially developed for AlphaLISA protein-protein interaction assays, in which the reaction requires at a low pH solution.
- Formulation:** AlphaLISA MES Buffer (5X) contains: 125 mM MES (pH=6.0), 1.25 M NaCl, , 0.5% Bovine serum albumin, 2.5% Triton X-100, 0.5 % Dextran and 0.25% Kathon CG II,
- Storage:** Store AlphaLISA MES Buffer (5X) at 4 °C.
- Reconstitution:** Add 4 volumes of Milli-Q™ grade water to 1 volume of 5X AlphaLISA MES Buffer.
- Stability:** The undiluted stock solutions are stable for at least 1 year from the manufacturing data when stored in original packing at 4 °C.

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

This product is not for resale or distribution except by authorized distributors.

LIMITED WARRANTY: PerkinElmer warrants that, at the time of shipment, the above named product is free from defects in material and workmanship and conforms to the specifications set forth above. PerkinElmer makes no other warranty, express or implied with respect to the product and expressly disclaims any warranty of merchantability or fitness for any particular purpose. Notification of any breach of the foregoing warranty must be made within 60 days of receipt of the product, unless otherwise provided in writing by PerkinElmer. No claim shall be honored if the customer fails to notify PerkinElmer within the period specified. The sole and exclusive remedy of the customer for any breach of the foregoing warranty is limited to either the replacement of the non-conforming product or the refund of the invoice price of the product. PERKINELMER SHALL NOT BE LIABLE FOR ANY DIRECT, INDIRECT, SPECIAL, INCIDENTAL, CONSEQUENTIAL OR PUNITIVE DAMAGES, WHETHER BASED ON CONTRACT, TORT, STRICT LIABILITY OR OTHERWISE, ARISING OUT OF THE DESIGN, MANUFACTURE, SALE, DELIVERY, OR USE OF THE PRODUCTS, EVEN IF THE LIMITED REMEDIES PROVIDED HEREIN FAIL OF THEIR ESSENTIAL PURPOSE OR PERKINELMER IS ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

PerkinElmer, Inc.
940 Winter Street
Waltham, MA 02451 USA
P: (800) 762-4000 or
(+1) 203-925-4602
www.perkinelmer.com

For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright© 2012, PerkinElmer, Inc. All rights reserved. PerkinElmer® is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners.