

AlphaLISA®

## Human KEAP1 Binding Kit

**Product number:** AL3194C                      **Lot Number:** 3255571

---

### Material provided:

**Kit Format:** AL3194C: 500 assay points                      AL3194F: 5000 assay points

The number of assay points is based on an assay volume of 20 µL in 384-well assay plates using kit components at the recommended concentrations.

**Standard Format:** 4.4 µg, lyophilized (one vial contains an amount of standard sufficient for performing 10 AlphaLISA standard curves)

**Manufacturing date:** 8/30/2023                      **Document version:** 1

---

### Product Information

**Application:** This kit is designed to select and characterize compounds that bind to human KEAP1 protein using a homogeneous AlphaLISA binding assay (no wash steps).

**Kit contents:** The kit contains 6 components: Anti-GST AlphaLISA Acceptor beads, Streptavidin-coated Donor beads, lyophilized Biotinylated KEAP1 ligand, lyophilized GST tagged KEAP1 protein, lyophilized KEAP1 standard (Ki696), and AlphaLISA Binding Assay Buffer.

**Storage:** Store kit in the dark at +4°C. Store reconstituted components at ≤-20°C at a minimum.

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

### Quality Control

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals and IC50 were measured on the EnVision Multilabel Plate Reader with Alpha option. We certify that these results meet our quality release criteria.

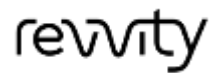
Maximum counts may vary between bead lots and the instrument used.

Maximum signal: 37042 counts

Minimum signal: 466 counts

IC50: 0.91 nM

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.

The Revvity logo consists of the word "revvity" in a lowercase, sans-serif font. The letters are black and have a slightly irregular, hand-drawn appearance.

**Revvity, Inc.**  
940 Winter Street  
Waltham, MA 02451 USA

(800) 762-4000 [www.revvity.com](http://www.revvity.com)

For a complete listing of our global offices, visit [www.revvity.com](http://www.revvity.com)  
Copyright ©2023, Revvity, Inc. All rights reserved.