

AlphaLISA®

KRAS G12C/SOS1 Binding Kit**Product number:** AL3151C**Lot Number:** 3275393**Material provided:****Kit Format:** AL3151C: 500 assay points AL3151F: 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.

Manufacturing date: 2/12/2024**Document version:** 1**Product Information**

Application: This kit is designed to assess inhibitors of KRAS G12C and SOS1 binding using a homogeneous no wash AlphaLISA assay.

Kit contents: The kit contains 6 components: anti-6xHis AlphaLISA Acceptor beads, Glutathione Alpha Donor beads, GTP, KRAS G12C protein, SOS1 protein and AlphaLISA PPI buffer.

Storage: The kit components must be stored at 4 °C in the dark. Reconstituted reagents can be aliquoted (not under 10µL) then frozen stored at – 80 °C for 1 month. Avoid multiple freeze-thaw cycles.

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

Quality Control

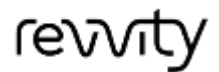
Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals and IC50 were measured on the EnVision Multilabel Plate Reader with Alpha option. We certify that these results meet our quality release criteria. Maximum counts may vary between bead lots and the instrument used.

IC50: 361.3 nM

SB: 106

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.

AL3151-R Rev01



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

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.