

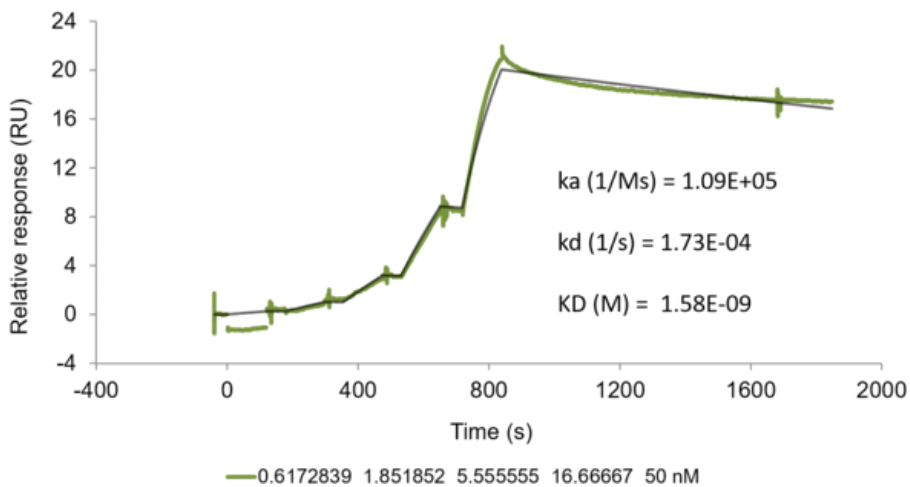
BCL-XL SPR Assay Service

Anti-apoptotic protein BCL-XL assay with [surface plasmon resonance \(SPR\)](#) technique

Target Overview

Target	Bcl-2-like protein 1 isoform L
HGNC symbol	BCL2L1
Synonyms	BCL-XL
Scientific information	Q07817 (https://www.uniprot.org/uniprot/Q07817)
Screening location	Malvern, PA, USA
Further information	More information can be found on our website SPR Assay Services for Drug Discovery

Binding kinetics of reference compound (ABT-737) to BCL-XL



The data was collected on a Biacore 8K using single-cycle kinetics. BCL-XL was immobilized to the sensor chip and the binding of ABT-737 was measured. The data was analyzed using a 1:1 kinetic binding model.

Let's Discover Together.

© May 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/bcl_xl_spr_malvern

