

## SPRINTer<sup>®</sup> K562 c-Myc Protein Turnover Biosensor Cell Line

**Catalog Number:** 91-1001C042

**Lot Number:** See Vial

**Contents:** 2 x 10<sup>6</sup> cells per vial in 1 mL

### Background

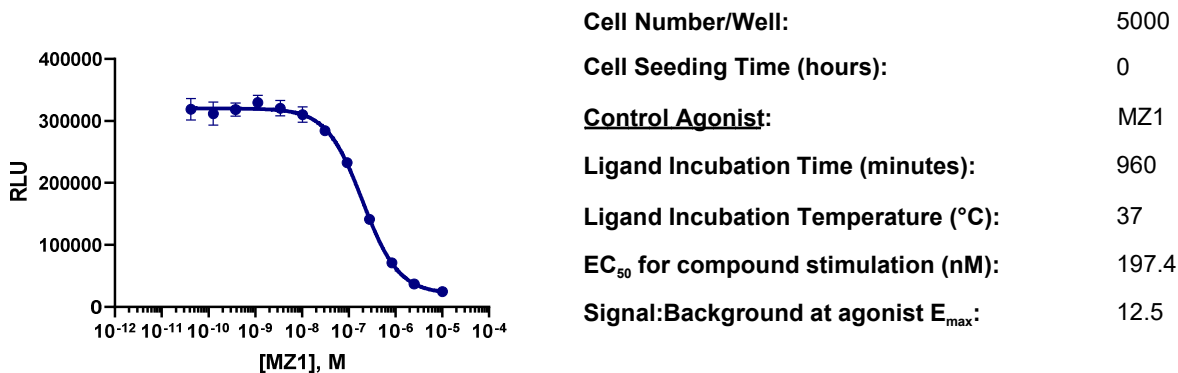
SPRINTer Protein Turnover Biosensor cell lines are engineered to introduce the EFC Enzyme Donor (ED) into the endogenous locus of the desired target gene. Expression of the target gene from its native promoter results in production of an ED-tagged target protein. Treatment of the engineered Biosensor cells with therapeutics that promote turnover of the target protein produces a decrease in EFC signal. Addition of exogenous Enzyme Acceptor (EA) and buffer lyses the cell and forces complementation of the ED and EA enzyme fragments. This results in the formation of a functional enzyme that hydrolyzes substrate to generate a chemiluminescent signal, allowing quantitation of drug-induced changes in ED-target protein levels.

### Product Information

**Cell Type:** K562  
**Species:** Human  
**Target Protein:** c-Myc  
**Description:** MYC proto-oncogene, bHLH transcription factor  
**Target Tag:** ePL  
**Tag Location:** N

### Functional Performance

Cells were plated in a 384-well plate and incubated at 37°C and 5% CO<sub>2</sub> to allow the cells to attach and grow. Cells were then stimulated with a control compound, using the assay conditions described below. Following stimulation, signal was detected using the PathHunter Detection Kit according to the recommended protocol. Please refer to page 2 for recommended assay reagents, detection reagents, and control compounds.



**Cell Number/Well:** 5000  
**Cell Seeding Time (hours):** 0  
**Control Agonist:** MZ1  
**Ligand Incubation Time (minutes):** 960  
**Ligand Incubation Temperature (°C):** 37  
**EC<sub>50</sub> for compound stimulation (nM):** 197.4  
**Signal:Background at agonist E<sub>max</sub>:** 12.5

**Recommended Culturing Conditions:** Start new cultures at 1 x 10<sup>5</sup> viable cells/mL; subculture when cells reach 1 x 10<sup>6</sup> cells/mL.

### Passage Stability

This cell line has been confirmed to be stable through a minimum of 10 passages with no significant drop in assay window or change in EC<sub>50</sub>.

### Mycoplasma Testing

This lot was tested and found to be free of mycoplasma contamination. Data available upon request.

### Required Materials

The following additional materials are required but not provided:

Product Use*	Product Description	Catalog Number
Detection	PathHunter® ProLabel®/ProLink™ Detection Kit	93-0812
Cell Culture	AssayComplete™ Cell Culture Kit-101	92-3101G
Cell Plating	AssayComplete™ Cell Plating 0 Reagent	93-0563R0A
Cell Detachment	AssayComplete™ Cell Detachment Reagent	92-0009
Cell Thawing	AssayComplete™ Thawing Reagent T6	92-4106TR
Cell Freezing	AssayComplete™ Freezing Reagent F5	92-5105FR

\*Please inquire about our cell line-specific AssayComplete Starter Packs to get you started with your cell culture needs.

### Required Antibiotics

Antibiotic Name	Concentration (µg/mL)	Catalog Number
AssayComplete™ Puromycin	Not Applicable	Not Applicable
AssayComplete™ Hygromycin B	Not Applicable	Not Applicable
AssayComplete™ G418	Not Applicable	Not Applicable

### Additional Ligand Information

**Control Agonist:** MZ1

**Vendor:** Tocris (Catalog No. 6154/5)

**Ordering:** +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
**Technical support:** +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
**General product information:** [www.discoverx.com](http://www.discoverx.com)

### Limited Use License Agreement

These products may be covered by issued US and/or foreign patents, patent application and subject to Limited Use Label License.

Please visit [discoverx.com/license](https://discoverx.com/license) for a list of products that are governed by limited use label license terms and relevant patent and trademark information.