
Optimization of the Tango™ RLN3R1-*bla* U2OS Cell Line

Tango™ RLN3R1-*bla* U2OS DA cells**Tango™ RLN3R1-*bla* U2OS cells**

Catalog Numbers – K1591 and K1569

Cell Line Descriptions

Tango™ RLN3R1-*bla* U2OS DA (Division Arrested) cells and Tango™ RLN3R1-*bla* U2OS cells contain the human Relaxin 3 receptor 1 (RLN3R1) linked to a TEV protease site and a Gal4-VP16 transcription factor stably integrated into the Tango™ GPCR-*bla* U2OS parental cell line. This parental cell line stably expresses a beta-arrestin/TEV protease fusion protein and the beta-lactamase (*bla*) reporter gene under the control of a UAS response element.

DA cells are irreversibly division arrested using a low-dose treatment of Mitomycin-C, and have no apparent toxicity or change in cellular signal transduction. Both the Tango™ RLN3R1-*bla* U2OS cells and the Tango™ RLN3R1-*bla* U2OS DA cells have been functionally validated for Z' factor and EC₅₀ concentrations of Relaxin-3 (Figure 1). In addition, Tango™ RLN3R1-*bla* U2OS cells have been tested for assay performance under variable conditions.

Validation Summary

Testing and validation of this assay was evaluated in a 384-well format using LiveBLAzer™-FRET B/G Substrate.

1. Relaxin-3 dose response under optimized conditions

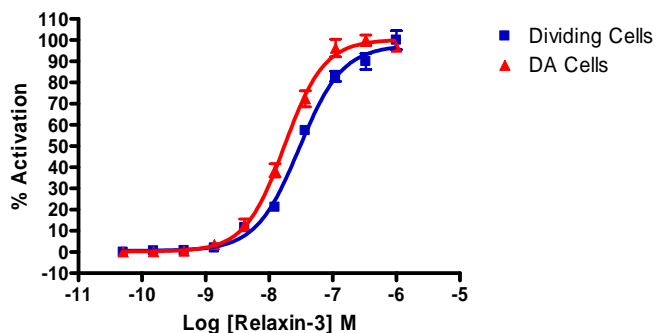
	<u>DA cells</u>	<u>Dividing Cells</u>
EC ₅₀	17.5nM	27.5nM
Z'-factor	0.85	0.73
Recommended cell no. /well	= 10,000	= 10,000
Recommended Stim. Time	= 5 hrs	= 5 hrs
Max. [Stimulation]	= 1000 nM	= 1000 nM

2. Antagonist dose response

No antagonists were commercially available at the time of publication of this document

Primary Agonist Dose Response

Figure 1 — Tango™ RLN3R1-bla U2OS cells and Tango™ RLN3R1-bla U2OS DA cells dose response to Relaxin-3 under optimized conditions



Tango™ RLN3R1-bla U2OS cells and Tango™ RLN3R1-bla U2OS DA cells (10,000 cells/well) were plated in a 384-well format and incubated for 16-20 hours. Cells were stimulated with a dilution series of Relaxin-3 (Phoenix Pharm. 035-36) in the presence of 0.1% DMSO for 5 hours. Cells were then loaded with LiveBLAzer™-FRET B/G Substrate for 2 hours. Fluorescence emission values at 460 nm and 530 nm were obtained using a standard fluorescence plate reader and % Activation plotted for each replicate against the concentrations of Relaxin-3.