
Optimization of the Tango™ HTR2A-*bla* U2OS Cell Line

Tango™ HTR2A-*bla* U2OS DA cells**Tango™ HTR2A-*bla* U2OS cells**

Catalog Numbers – K1613 and K1562

Cell Line Descriptions

Tango™ HTR2A-*bla* U2OS DA (Division Arrested) cells and Tango™ HTR2A-*bla* U2OS cells contain the human Serotonin Type 2a receptor (HTR2A) 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™ HTR2A-*bla* U2OS cells and the Tango™ HTR2A-*bla* U2OS DA cells have been functionally validated for Z' factor and EC₅₀ concentrations of 5-Hydroxytryptamine (Figure 1). In addition, Tango™ HTR2A-*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. 5-Hydroxytryptamine dose response under optimized conditions

	DA cells	Dividing Cells
EC ₅₀	16.29 nM	17.9 nM
Z'-factor	0.72	0.85
Recommended cell no. /well		= 10,000
Recommended Stim. Time		= 5 hrs
Max. [Stimulation]		= 10000 nM

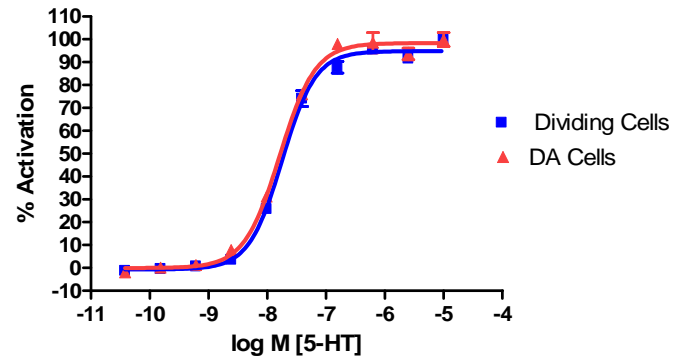
2. Antagonist dose response

Mianserin:

Dividing IC ₅₀	= 28 nM
Cryopreserved IC ₅₀	= 68 nM
Division Arrested IC ₅₀	= 58 nM

Primary Agonist Dose Response

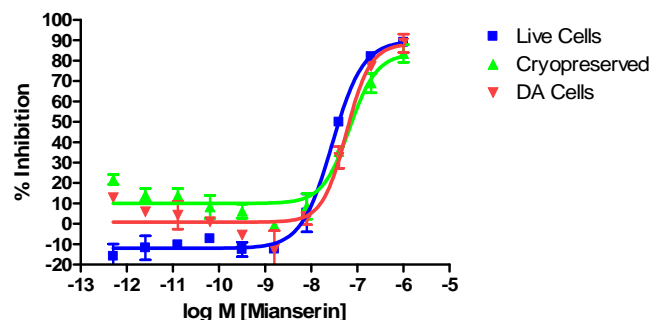
Figure 1 — Tango™ HTR2A-*bla* U2OS cells and Tango™ HTR2A-*bla* U2OS DA cells dose response to 5-Hydroxytryptamine under optimized conditions



Tango™ HTR2A-*bla* U2OS cells and Tango™ HTR2A-*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 5-Hydroxytryptamine (Sigma H9523) 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 5-Hydroxytryptamine.

Antagonist Dose Response

Figure 2 — Tango™ HTR2A-*bla* U2OS cells dose response to Mianserin



Tango™ HTR2A-*bla* U2OS cells (10,000 cells/well) were plated in a 384-well format and incubated for 16-20 hours. Cells were exposed to Mianserin (Sigma M2525) for 30 min. and then stimulated with an EC80 concentration of 5-Hydroxytryptamine (Sigma H9523) 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 for the various substrate loading times were obtained using a standard fluorescence plate reader and the % Inhibition plotted against the indicated concentrations of Mianserin.