
Optimization of the Tango™ MTNR1A-*bla* U2OS Cell Line

Tango™ MTNR1A-*bla* U2OS DA Assay Kit**Tango™ MTNR1A-*bla* U2OS cells**

Catalog Numbers – K1670 and K1669

Cell Line Descriptions

Tango™ MTNR1A-*bla* U2OS DA (Division Arrested) cells and Tango™ MTNR1A-*bla* U2OS cells contain the human Melatonin Receptor 1A (MTNR1A) 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 reporter gene under the control of a UAS response element. Division Arrested (DA) cells are available in an Assay Kit, which includes cells and sufficient substrate to analyze 1 x 384-well plate.

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™ MTNR1A-*bla* U2OS cells and the Tango™ MTNR1A-*bla* U2OS DA cells have been functionally validated for Z' factor and EC₅₀ concentrations of Melatonin (Figure 1). In addition, Tango™ MTNR1A-*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. Melatonin dose response under optimized conditions

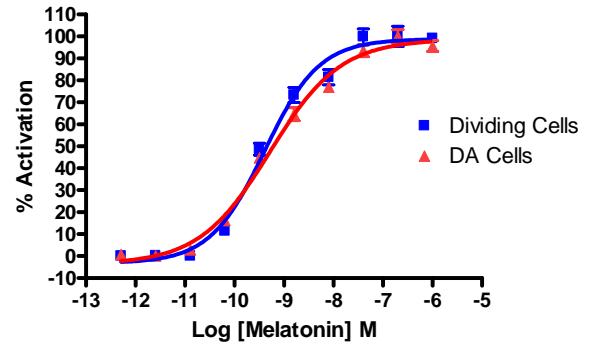
| | DA cells | Dividing Cells |
|----------------------------|-----------|----------------|
| EC ₅₀ | 519 pM | 404 pM |
| Z'-factor | 0.78 | 0.77 |
| Recommended cell no. /well | = 10,000 | = 10,000 |
| Recommended Stim. Time | = 5 hrs | = 5 hrs |
| Max. [Stimulation] | = 1000 nM | = 1000 nM |

2. Antagonist dose response

Luzindole IC₅₀ = 702 nM

Primary Agonist Dose Response

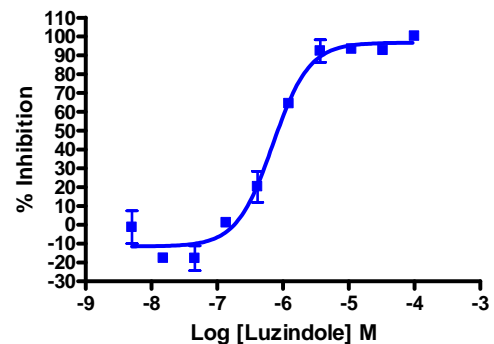
Figure 1 — Tango™ MTNR1A-*bla* U2OS cells and Tango™ MTNR1A-*bla* U2OS DA cells dose response to Melatonin under optimized conditions



Tango™ MTNR1A-*bla* U2OS cells and Tango™ MTNR1A-*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 Melatonin (Sigma M5250) 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 Melatonin.

Antagonist Dose Response

Figure 3 — Tango™ MTNR1A-*bla* U2OS cells dose response to Luzindole



Tango™ MTNR1A-*bla* U2OS cells (10,000 cells/well) were plated in a 384-well format and incubated for 16-20 hours. Cells were exposed to Luzindole (Sigma L2407) for 30 min. and then stimulated with an EC₈₀ concentration of Melatonin (Sigma M5250) 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 Luzindole.