
Optimization of the Tango™ D1-*bla* U2OS Cell Line

Tango™ D1-*bla* U2OS DA cells**Tango™ D1-*bla* U2OS cells**

Catalog Numbers – K1505 and K1485

Cell Line Descriptions

Tango™ D1-*bla* U2OS DA (Division Arrested) cells and Tango™ D1-*bla* U2OS cells contain the human Dopamine Receptor 1 (D1) 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™ D1-*bla* U2OS cells and the Tango™ D1-*bla* U2OS DA cells have been functionally validated for Z' factor and EC₅₀ concentrations of Dihydropyridine (Figure 1). In addition, Tango™ D1-*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. Dihydraxidine dose response under optimized conditions

| | DA cells | Dividing Cells |
|----------------------------|-------------|----------------|
| EC ₅₀ | 400 nM | 249 nM |
| Z'-factor | 0.77 | 0.76 |
| Recommended cell no. /well | = 15,000 | = 15,000 |
| Recommended Stim. Time | = 5 hrs | = 5 hrs |
| Max. [Stimulation] | = 25,000 nM | = 25,000 nM |

2. Antagonist dose response

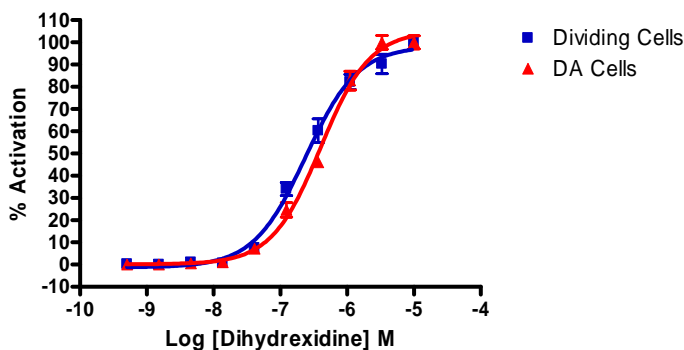
R(+)-SCH-23390 IC₅₀ = 2.1 nM

3. Agonist 2nd messenger dose response.

Dihydraxidine EC₅₀ = 15 nM

Primary Agonist Dose Response

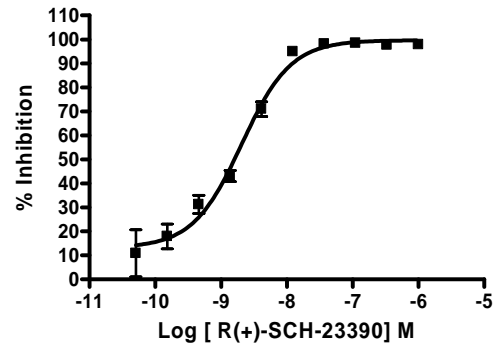
Figure 1 — Tango™ D1-*bla* U2OS cells and Tango™ D1-*bla* U2OS DA cells dose response to Dihydraxidine under optimized conditions



Tango™ D1-*bla* U2OS cells and Tango™ D1-*bla* U2OS DA cells (15,000 cells/well) were plated in a 384-well format and incubated for 16-20 hours. Cells were stimulated with a dilution series of Dihydraxidine (Tocris 884) 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.

Antagonist Dose Response

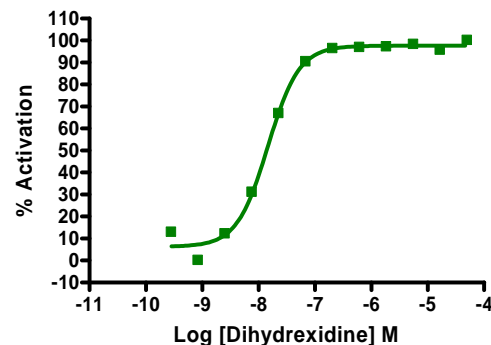
Figure 2 — Tango™ D1-*bla* U2OS cells dose response to R (+)-SCH-23390



Tango™ D1-*bla* U2OS cells (15,000 cells/well) were plated in a 384-well format and incubated for 16-20 hours. Cells were exposed to () for 30 min. and then stimulated with an EC80 concentration of Dihydraxidine (Tocris 884) 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 R(+)-SCH-23390

2nd Messenger Dose Response

Figure 3 — Tango D1-*bla* U2OS cells 2nd messenger dose response to Dihydraxidine under optimized conditions.



Tango™ D1-*bla* U2OS cells were tested for a response to dihydraxidine with a TR-FRET cAMP kit.