

---

**Optimization of the Tango™ CCR5-*bla* U2OS Cell Line**

---

**Tango™ CCR5-*bla* U2OS cells**

Catalog Numbers – K1788

**Cell Line Descriptions**

The Tango™ CCR5-*bla* U2OS cells contain the human Chemokine (C-C Motif) Receptor 5 (CCR5) 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.

The Tango™ CCR5-*bla* U2OS cells have been functionally validated for Z' factor and EC<sub>50</sub> concentrations of MIP-1a (Figure 1). In addition, Tango™ CCR5-*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. MIP-1a dose response under optimized conditions

|                            |                       |
|----------------------------|-----------------------|
|                            | <u>Dividing Cells</u> |
| EC <sub>50</sub>           | 18.5 nM               |
| Z'-factor                  | 0.89                  |
| Recommended cell no. /well | = 10,000              |
| Recommended Stim. Time     | = 16 hrs              |
| Max. [Stimulation]         | = 641 nM              |

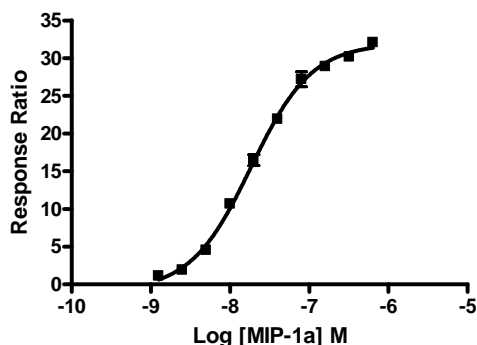
### 2. Assay performance with variable cell number.

### 3. Antagonist dose response

IC<sub>50</sub> = 324 nM

## Primary Agonist Dose Response

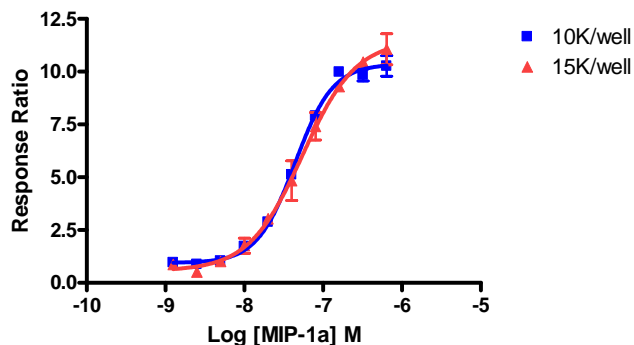
Figure 1 — Tango™ CCR5-bla U2OS cells dose response to MIP-1a under optimized conditions



Tango™ CCR5-bla U2OS 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 MIP-1a (Biosource (IVGN) PHC1105) in the presence of 0.1% DMSO for 16 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 Response Ratios plotted for each replicate against the concentrations of MIP-1a.

## Assay Performance with Variable Cell Number

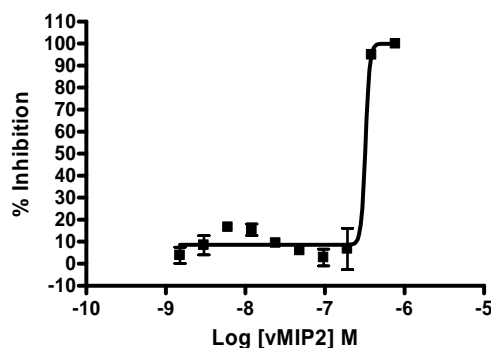
Figure 2 — Tango™ CCR5-bla U2OS cells dose response to MIP-1a with 10K or 15K cells/well



Tango™ CCR5-bla U2OS cells were plated in a 384-well format at 10,000 or 15,000 cells/well and incubated for 16-24 hours. On the day of the assay, cells were stimulated with MIP-1a (Biosource (IVGN) PHC1105) in the presence of 0.1% DMSO for 16 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 Response Ratios plotted for each replicate against the concentrations of MIP-1a.

## Antagonist Dose Response

Figure 3 — Tango™ CCR5-bla U2OS cells dose response to vMIP2



Tango™ CCR5-bla U2OS cells (10,000 cells/well) were plated in a 384-well format and incubated for 16-20 hours. Cells were exposed to vMIP2 for 30 min. and then stimulated with an EC80 concentration of MIP-1a (Biosource (IVGN) PHC1105) in the presence of 0.1% DMSO for 16 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 not currently available.