
Optimization of the Tango™ CXCR6-*bla* U2OS Cell Line

Tango™ CXCR6-*bla* U2OS DA cells**Tango™ CXCR6-*bla* U2OS cells**

Catalog Numbers – K1583 and K1563

Cell Line Descriptions

Tango™ CXCR6-*bla* U2OS DA (Division Arrested) cells and Tango™ CXCR6-*bla* U2OS cells contain the human Chemokine (C-X-C motif) receptor 6 (CXCR6) 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™ CXCR6-*bla* U2OS cells and the Tango™ CXCR6-*bla* U2OS DA cells have been functionally validated for Z' factor and EC₅₀ concentrations of CXCL16 (Figure 1).

Validation Summary

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

1. CXCL16 dose response under optimized conditions

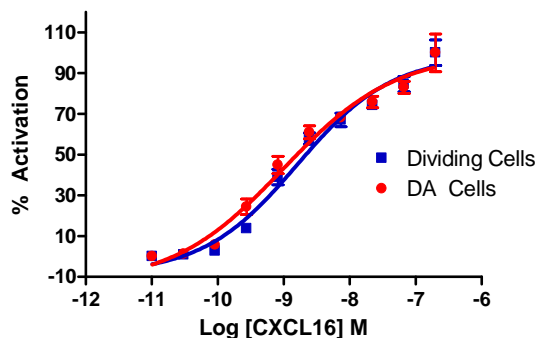
	DA cells	Dividing Cells
EC ₅₀	1.0 nM	1.6 nM
Z'-factor	0.71	0.66
Recommended cell no. /well	= 10,000	= 10,000
Recommended Stim. Time	= 5 hrs	= 5 hrs
Max. [Stimulation]	= 200 nM	= 200 nM

2. Antagonist dose response

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

Primary Agonist Dose Response

Figure 1 — Tango™ CXCR6-bla U2OS cells and Tango™ CXCR6-bla U2OS DA cells dose response to CXCL16 under optimized conditions



Tango™ CXCR6-bla U2OS cells and Tango™ CXCR6-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 CXCL16 (Sigma C8615) 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 plotted for each replicate against the concentrations of CXCL16.