

## MULTISCREEN™ DIVISION ARRESTED CELL LINE HEK293T $\beta$ -Arrestin1

### Data sheet

#### PRODUCT INFORMATION

**Catalog Number:** DCA0001BA1

**Lot Number:** DCA0001BA1-031326

**Quantity:** 1 vial ( $4 \times 10^6$ ) frozen cells

**Freeze Medium:** Cellbanker 2

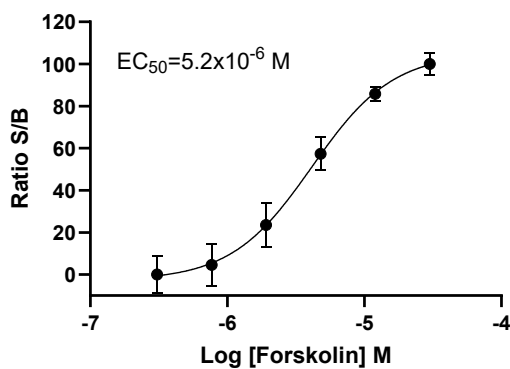
**Host cell:** HEK293T  $\beta$ -Arrestin1

**Recommended Storage:** Liquid nitrogen upon receiving

**Propagation Medium:** DMEM, 10% FBS

**Application:** Functional assays

**Figure 1**



**Figure 1.** Dose-dependent stimulation of intracellular cAMP accumulation by forskolin, measured with MULTISCREEN™ TR-FRET cAMP 1.0 No Wash Assay Kit (Multispan MSCM01).

**FOR RESEARCH USE ONLY.**

All rights reserved. No part of this document may be reproduced in any form without prior permission in writing