



Cell transient transfection with tagged GPCR or RTK plasmids

Revvity proposes a broad list of GPCR and RTK plasmids encoding SNAP-tag®, CLIP-tag® and HaloTag® at the N-terminal position. Once transfected, cells expressing the tagged receptor can be labeled with the specific substrate derivatized with HTRF fluorophores.

Material

Plasmids

- The list of GPCR and TRK plasmids (human) is available on our website.

Recommended reagents

- HEK 293 cells
- PBS (Life technologies Cat.# 10010-031)
- Cell dissociation buffer (Sigma Cat.#C5914)
- Opti-MEM, with GlutaMAX and phenol red (Life technologies Cat.# 51985-026)
- Lipofectamine 2000 (Life technologies Cat.#11668-027)
- Culture medium for HEK293: DMEM (Life technologies Cat.#31966-021) supplemented with Penicillin-Streptomycin 1X (Life technologies Cat.# 15070-063), HEPES 2mM (Life technologies Cat.#15630-122), MEM 1X (Life Technologies Cat.#11140-035)

Transient transfection - flask T175

Two to three days before the transfection, seed the cells in a T175 flask. For an efficient transfection, cells should be adherent with a confluency around 70 - 80%

1. Prepare the transfection mix, add in the following order

- Opti-NEM: 8 mL
- Lipofectamine: 60 µL
- Plasmid (1µg/µL): 20 µL

Incubate for 20 minutes at RT

2. Gently remove the cell culture medium by aspiration
3. Wash cells with 5 mL of PBS
4. Add the 8.08 mL of transfection mix preparation then add 12 mL of cell culture medium
5. Incubate for 1 night at 37°C + 5% CO₂.
6. Cells expressing the substrate-tag receptor can now be labelled with Tag-lite SNAP-Lumi4-Tb and used to carry out a binding assay



The information provided in this document is for reference purposes only and may not be all-inclusive. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume liability for the accuracy or completeness of the information contained herein. Users should exercise caution when handling materials as they may present unknown hazards. Revvity shall not be liable for any damages or losses resulting from handling or contact with the product, as Revvity cannot control actual methods, volumes, or conditions of use. Users are responsible for ensuring the product's suitability for their specific application. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF

Manufactured by Cisbio Bioassays - Parc Marcel Boiteux - 30200 Codolet - FRANCE

www.revvity.com

revvity