

## CELLULAR FUNCTIONAL GPCR ASSAYS

Functional GPCR assays are cell-based assays to measure agonist-like and antagonist activity of compounds. Various technologies are used: TR-FRET to determine cAMP concentration and IP1 levels, real time fluorescence to monitor calcium flux, cellular dielectric spectroscopy to measure impedance modulation. Large batches of cells, from which whole cells are frozen and stored for functional assays, are produced in-house. Cells are quality controlled before use. A novel patented host cell line was designed and constructed for Gi protein coupled receptors.

### NON-PEPTIDE RECEPTORS

Family/assay	Ref.	Family/assay	Ref.
<b>ADENOSINE</b>		<b>LYSOPHOSPHOLIPID</b>	
A <sub>1</sub>	G127	Lysophosphatidic acid	
A <sub>2A</sub>	G002	LPA <sub>2</sub>	G144
A <sub>2B</sub>	G003	LPA <sub>3</sub>	G145
A <sub>3</sub>	G107	Sphingosine 1-phosphate	
<b>ADRENERGIC</b>		S <sub>1</sub> P <sub>2</sub>	G026
alpha <sub>1A</sub>	G004	S <sub>1</sub> P <sub>3</sub>	G027
alpha <sub>1B</sub>	G005	<b>MELATONIN</b>	
alpha <sub>2A</sub>	G120	MT <sub>1</sub> (ML <sub>1A</sub> )	G136
alpha <sub>2B</sub>	G011	MT <sub>2</sub> (ML <sub>1B</sub> )	G028
alpha <sub>2C</sub>	G006	<b>MUSCARINIC</b>	
beta <sub>1</sub>	G007	M <sub>1</sub>	G029
beta <sub>2</sub>	G008	M <sub>2</sub>	G030
beta <sub>3</sub>	G009	M <sub>3</sub>	G031
<b>CALCIUM SENSING</b>		M <sub>4</sub>	G032
CaS	G010	M <sub>5</sub>	G033
<b>CANNABINOID</b>		<b>PLATELET ACTIVATING FACTOR</b>	
CB <sub>1</sub>	G012	PAF	G034
CB <sub>2</sub>	G013	<b>PROSTANOID</b>	
<b>DOPAMINE</b>		DP	G121
D <sub>1</sub>	G014	EP <sub>1</sub>	G035
D <sub>2S</sub>	G116	EP <sub>2</sub>	G036
D <sub>3</sub>	G016	EP <sub>3</sub>	G122
D <sub>4.4</sub>	G017	EP <sub>4</sub>	G037
D <sub>5</sub>	G018	FP	G038
<b>FREE FATTY ACID</b>		IP (PGI <sub>2</sub> )	G039
FFA1 (GPR40)	G123	TP (TXA <sub>2</sub> /PGH <sub>2</sub> )	G040
FFA2 (GPR43)	G132	<b>PURINERGIC</b>	
FFA3 (GPR41)	G143	P2Y <sub>1</sub>	G041
<b>GABA</b>		P2Y <sub>2</sub>	G108
GABA <sub>B</sub>	G111	P2Y <sub>4</sub>	G042
<b>HISTAMINE</b>		P2Y <sub>6</sub>	G043
H <sub>1</sub>	G115	<b>SEROTONIN</b>	
H <sub>2</sub>	G020	5-HT <sub>1A</sub>	G044
H <sub>3</sub>	G021	5-HT <sub>1B</sub>	G126
H <sub>4</sub>	G125	5-HT <sub>1D</sub>	G112
<b>LEUKOTRIENES</b>		5-HT <sub>2A</sub>	G046
BLT <sub>1</sub> (LTB <sub>4</sub> )	G023	5-HT <sub>2B</sub>	G047
CysLT <sub>1</sub> (LTD <sub>4</sub> )	G024	5-HT <sub>2C</sub>	G048
CysLT <sub>2</sub> (LTC <sub>4</sub> )	G025	5-HT <sub>4e</sub>	G049
		5-HT <sub>6</sub>	G050
		5-HT <sub>7</sub>	G051

## PEPTIDE RECEPTORS

Family/assay	Ref.
<b>ANGIOTENSIN II</b>	
☐ AT <sub>1</sub>	🇫 G052
<b>APELIN</b>	
☐ APJ (apelin)	🇫 G146
<b>BOMBESIN</b>	
☐ BB <sub>1</sub>	🇫 G147
☐ BB <sub>2</sub>	🇫 G053
☐ BB <sub>3</sub>	🇫 G054
<b>BRADYKININ</b>	
☐ B <sub>1</sub>	🇫 G055
☐ B <sub>2</sub>	🇫 G056
<b>CALCITONIN</b>	
☐ CT (calcitonin)	🇫 G057
<b>CALCITONIN GENE-RELATED PEPTIDE</b>	
☐ CGRP	🇫 G058
<b>CHEMOKINES</b>	
☐ CCR1	🇫 G059
☐ CCR2	🇫 G138
☐ CCR3	🇫 G148
☐ CXCR2 (IL-8B)	🇫 G113
☐ CXCR4	🇫 G114
<b>CHOLECYSTOKININ</b>	
☐ CCK <sub>1</sub> (CCK <sub>A</sub> )	🇫 G062
☐ CCK <sub>2</sub> (CCK <sub>B</sub> )	🇫 G063
<b>COMPLEMENT 5A</b>	
☐ C5α	🇫 G064
<b>CORTICOTROPIN RELEASING FACTOR</b>	
☐ CRF <sub>1</sub>	🇫 G065
☐ CRF <sub>2α</sub>	🇫 G066
<b>ENDOTHELIN</b>	
☐ ET <sub>A</sub>	🇫 G067
☐ ET <sub>B</sub>	🇫 G068
<b>N-FORMYL PEPTIDE</b>	
☐ fMLP	🇫 G069
<b>GALANIN</b>	
☐ GAL <sub>2</sub>	🇫 G070
<b>GLUCAGON</b>	
☐ GLP-1	G072
☐ GLP-2	🇫 G073
☐ glucagon	🇫 G074
☐ secretin	🇫 G075
<b>GLYCOPROTEIN HORMONE</b>	
☐ TSH	🇫 G076
<b>GROWTH HORMONE-RELEASING HORMONE</b>	
☐ GHRH	🇫 G071
<b>MELANIN CONCENTRATING HORMONE</b>	
☐ MCH <sub>1</sub>	🇫 G077
☐ MCH <sub>2</sub>	🇫 G078

Family/assay	Ref.
<b>MELANOCORTIN</b>	
☐ MC <sub>1</sub>	G079
☐ MC <sub>2</sub>	🇫 G080
☐ MC <sub>3</sub>	🇫 G081
☐ MC <sub>4</sub>	🇫 G082
☐ MC <sub>5</sub>	🇫 G083
<b>MOTILIN</b>	
☐ motilin	🇫 G084
<b>NEUROKININ</b>	
☐ NK <sub>1</sub>	🇫 G085
☐ NK <sub>2</sub>	🇫 G086
☐ NK <sub>3</sub>	🇫 G087
<b>NEUROPEPTIDE Y</b>	
☐ Y <sub>1</sub>	🇫 G088
<b>NEUROTENSIN</b>	
☐ NTS <sub>1</sub> (NT <sub>1</sub> )	🇫 G089
<b>OPIOID AND OPIOID-LIKE</b>	
☐ delta <sub>2</sub> (DOP)	G129
☐ kappa (KOP)	G090
☐ mu (MOP)	🇫 G091
☐ NOP (ORL1)	🇫 G119
<b>OREXIN</b>	
☐ OX <sub>1</sub>	🇫 G092
☐ OX <sub>2</sub>	🇫 G093
<b>PARATHYROID HORMONE</b>	
☐ PTH1	🇫 G130
<b>PROKINETICIN</b>	
☐ PK <sub>1</sub>	🇫 G117
☐ PK <sub>2</sub>	🇫 G118
<b>PROTEINASE-ACTIVATED</b>	
☐ PAR1	🇫 G109
☐ PAR2	🇫 G110
<b>RELAXIN</b>	
☐ RXFP1	🇫 G094
<b>SOMATOSTATIN</b>	
☐ sst <sub>1</sub>	🇫 G095
☐ sst <sub>4</sub>	🇫 G096
☐ sst <sub>5</sub>	🇫 G097
<b>THYROTROPIN RELEASING HORMONE</b>	
☐ TRH <sub>1</sub>	🇫 G098
<b>UROTENSIN-II</b>	
☐ UT	🇫 G099
<b>VASOACTIVE INTESTINAL PEPTIDE</b>	
☐ PAC <sub>1</sub> (PACAP)	🇫 G100
☐ VPAC <sub>1</sub> (VIP <sub>1</sub> )	🇫 G101
☐ VPAC <sub>2</sub> (VIP <sub>2</sub> )	🇫 G102
<b>VASOPRESSIN</b>	
☐ OT	🇫 G103
☐ V <sub>1α</sub>	🇫 G104
☐ V <sub>2</sub>	🇫 G105

☐ cellular assay ☐ new assay ☐ new protocol 🇫 human

## TESTING CONDITIONS

### SUGGESTED TESTING

Primary screening on both agonist and antagonist effects at 1-10  $\mu\text{M}$  in duplicate (2 wells), followed-up for  $\text{IC}_{50}/\text{K}_{\text{Bapp}}$  or  $\text{EC}_{50}$  determination (8 concentrations in duplicate (16 wells) when compound displays more than 50% inhibition or 50% of agonist response.

### SAMPLE SIZE (including $\text{IC}_{50}$ follow-up studies)

Assuming a molecular weight  $\leq 500$  g/mol and a testing concentration of 10  $\mu\text{M}$  in duplicate (including a possible retest).

	SCREENING		SCREENING + FOLLOW UP <sup>1</sup>	
	WEIGHT (pre-weighed)	VOLUME (100% DMSO)	WEIGHT (pre-weighed)	VOLUME (100% DMSO)
<b>INDIVIDUAL CATALOG ASSAYS</b>				
1 to 3 assays	1 mg	25 $\mu\text{L}$ @ 10 mM	1 mg	60 $\mu\text{L}$ @ 10 mM
4 to 5 assays	1 mg	35 $\mu\text{L}$ @ 10 mM	1 mg	70 $\mu\text{L}$ @ 10 mM
6 to 10 assays	1 mg	50 $\mu\text{L}$ @ 10 mM	1 mg	100 $\mu\text{L}$ @ 10 mM
11 to 15 assays	1 mg	65 $\mu\text{L}$ @ 10 mM	1 mg	110 $\mu\text{L}$ @ 10 mM
16 to 20 assays	1 mg	75 $\mu\text{L}$ @ 10 mM	1 mg	125 $\mu\text{L}$ @ 10 mM
21 to 40 assays	1 mg	100 $\mu\text{L}$ @ 10 mM	1.5 mg	250 $\mu\text{L}$ @ 10 mM
41 to 50 assays	1 mg	150 $\mu\text{L}$ @ 10 mM	2 mg	275 $\mu\text{L}$ @ 10 mM
51 to 70 assays	1.5 mg	225 $\mu\text{L}$ @ 10 mM	2.5 mg	400 $\mu\text{L}$ @ 10 mM
71 to 100 assays	2 mg	300 $\mu\text{L}$ @ 10 mM	3 mg	550 $\mu\text{L}$ @ 10 mM
101 to 135 assays	2 mg	350 $\mu\text{L}$ @ 10 mM	4 mg	650 $\mu\text{L}$ @ 10 mM
136 to 150 assays	2.5 mg	400 $\mu\text{L}$ @ 10 mM	4 mg	750 $\mu\text{L}$ @ 10 mM
151 to 200 assays	3 mg	500 $\mu\text{L}$ @ 10 mM	5 mg	1000 $\mu\text{L}$ @ 10 mM
201 to 250 assays	3.5 mg	600 $\mu\text{L}$ @ 10 mM	inquire	inquire

<sup>1</sup> Assuming ~10% of test in  $\text{EC}_{50}$  or  $\text{IC}_{50}$ . Usually, for 1 curve: 30  $\mu\text{L}$ @ 10 mM and + 25  $\mu\text{L}$ @ 10 mM by additional  $\text{EC}_{50}$  or  $\text{IC}_{50}$  determination.

### REQUESTED COMPOUND INFORMATION

To reduce the registration time and ensure that all the appropriate information is available to start the study in a shortest possible timeframe, please use **Cerep compound submission form** <sup>1</sup> or MS Excel file <sup>2</sup>, and provide the following **compound information**:

- **Name (compound ID) / Batch # / Molecular weight** <sup>3</sup> / **Formula weight** <sup>4</sup> / **Stock concentration** / **Stock solvent** / **Quantity** / **Unit** <sup>5</sup> / **Form** / **Storage conditions** / **Solubility**, as well as **Plate ID/plate position** for compounds delivered in plates, **Comments** <sup>6</sup>, and **Quotation number**.

**NOTE:** Impurity and colored compounds might affect the results (compound color information is mentioned in the study report).

#### ► General remarks:

- **If compound(s) are supplied as a stock solution in plate(s)** (preferred format for any submission of 10 or more compounds), please leave columns 1 and 12 empty in a 96W plate. The 384W plate format is also acceptable with columns 1, 2, 23 and 24 empty. For any other plate format, please inquire.
- **If compound(s) are not soluble in 100% DMSO**, please provide any useful information concerning the solubility of the compound. The following solvents are compatible with most of our assays: DMSO (Cerep standard),  $\text{H}_2\text{O}$ , Methanol, Tris/HCl 10 mM pH 7.4.
- **Organic solvents such as acetone, chloroform, ether, acetonitrile, tetrahydrofuran and trifluoroacetic acid are not recommended** as they will significantly affect the results from many *in vitro* assays, even at very low concentrations.

**WARNING:** Cerep will apply the standard solubilization process when compounds are received at the testing site, unless special instructions are provided with the compounds.

Customized handling procedure of compounds can be accommodated, please inquire for pricing conditions.

### PROTOCOL

A **typical protocol for all cellular assays** includes a minimum of 6-control wells (background, and maximal signal with and without vehicle) plus an 8-point dose-response of the relevant reference compound.

The reference compound for each assay is listed in each assay description. The historical average  $\text{EC}_{50}$  or  $\text{IC}_{50}$  value is also shown in each assay description.

► ANY OF OUR ASSAY PROTOCOLS CAN BE CUSTOMIZED: PLEASE INQUIRE

<sup>1</sup> Cerep compound submission form will be emailed to you with your quotation. A copy can be requested from sales@cerep.com, or downloaded from Cerep website: www.cerep.com/Catalog Online

<sup>2</sup> Systematically required for studies of 10 compounds or more.

<sup>3</sup> Molecular weight (MW) of free acid or base form.

<sup>4</sup> Formula weight (FW) including salt form and/or hydrate form if applicable.

<sup>5</sup> mg $\text{g}^{-1}$ , mL $\text{g}^{-1}$

<sup>6</sup> e.g. useful information such as sensitivity to light, stability or hygroscopicity issues.

## ■ DELIVERABLES

Percent effect or percent inhibition (mean of replicates), individual values as percent of control, EC<sub>50</sub> or IC<sub>50</sub> value (calculated from a minimum 5 concentration testing), K<sub>Bapp</sub>, Hill coefficient (nH), and plotted EC<sub>50</sub> or IC<sub>50</sub> curves.

## ■ DATA TURNAROUND

**Complete data set is typically available within 3 weeks**, after receipt of the compounds at the testing site (providing that we receive all available information to initiate the study).

Secure, password-protected data can be viewed on line as soon as they are produced, after scientific approval and QC-ed by an experienced technician.

► FOR MAJORITY OF ASSAYS, EXPRESSCREEN OPTION (5 BUSINESS DAYS) IS AVAILABLE:  
PLEASE INQUIRE



### FRANCE

Le Bois l'Evêque  
86600 CELLE L'EVESCAULT  
tel. +33 (0)5 49 89 30 00

(Headquarters)

155 boulevard Haussmann  
75008 PARIS  
tel. +33 (0)1 45 64 44 60

### USA

15318 N.E. 95th Street  
REDMOND, WA 98052  
tel. +1 (425) 895 8666

### JAPAN

Namiki Shoji Co., Ltd.  
Kenseishinjuku Bldg. 5-5-3  
Shinjuku, Shinjuku-ku  
TOKYO, 160-0022  
tel. +81 (0)3 3354 4026  
fax +81 (0)3 3352 2196

### CHINA

Ai Di Sheng (Edison) Road 326,  
302-1 room  
Zhangjiang High-Tech Park  
SHANGHAI  
tel. +86 18702160370

## ASSAY CATALOG REFERENCES

In 2009, Cerep finalized one of the milestones in the industrialization process: the implementation of new referencing and supply chain management systems. The references of each of the assays have thus been simplified: **each reference will now be displayed as 4 digits.**

A correlation table between old and new assay references is available at [www.cerep.com](http://www.cerep.com) **CATALOG ONLINE**  
<<http://www.cerep.com/Cerep/Users/pages/Catalog/Assay/catalog.asp>>

## QUESTIONS OR CONCERNS?

Please contact us: [sales@cerep.com](mailto:sales@cerep.com)