




TISSUE ASSAYS

Tissue bioassays are functional assays designed to evaluate the agonist and antagonist activities of compounds at various receptors and ion channels in whole tissues. The tissues used are isolated from contractile cardiac and smooth muscles of the cardiovascular, respiratory, gastrointestinal and urogenital tracts of rodents, rabbit or human. The selected tissues are assayed in organ bath systems to record the contractile activity in conditions as selective as possible to avoid any potential interaction with other receptors known to be present in the tissue used.

NON-PEPTIDE RECEPTORS ■

Family/assay	Ref.
ADENOSINE	
T A ₁	0294
T A _{2A}	0295
T A _{2B}	0783
ADRENERGIC	
T alpha ₁ (non-selective)	0296
T alpha ₂ (non-selective)	0297
T beta ₁	0298 
T beta ₂	0299
T beta ₃	0300
CANNABINOID	
T CB ₁	0307
DOPAMINE	
T D ₁	0309
T D ₂	0310
GABA	
T GABA _A	0314
T GABA _B	0315
HISTAMINE	
T H ₁	0316
T H ₂	0317
MUSCARINIC	
T M ₁	0318
T M ₂	0319
T M ₃	0320
PURINERGIC	
T P2X	0332
SEROTONIN	
T 5-HT _{2A}	0334
T 5-HT _{2B}	0333
T 5-HT ₃	0335
T 5-HT ₄	0336
SIGMA	
T sigma (non-selective)	0337

PEPTIDE RECEPTORS ■

Family/assay	Ref.
ANGIOTENSIN II	
T AT ₁	0301
BOMBESIN	
T BB ₁	0303
BRADYKININ	
T B ₁	0935 
T B ₁	0304
T B ₂	0623 
T B ₂	0305
CALCITONIN GENE-RELATED PEPTIDE	
T CGRP	0306
CHOLECYSTOKININ	
T CCK ₁ (CCK _A)	0308

Family/assay	Ref.
ENDOTHELIN	
T ET _A	0311
T ET _B	0312
NEUROKININ	
T NK ₁	0321
T NK ₁	0624
T NK ₂	0322
T NK ₃	0323
T NK ₃	0625
NEUROPEPTIDE Y	
T Y ₁	0324
T Y ₂	0325
T Y ₃	0326
NEUROTENSIN	
T NT (non-selective)	0327
OPIOID AND OPIOID-LIKE	
T delta ₂ (DOP)	0328
T kappa (KOP)	0329
T mu (MOP)	0330
T NOP (ORL1)	1226
VASOACTIVE INTESTINAL PEPTIDE	
T VPAC ₂ (VIP ₂)	0622
VASOPRESSIN	
T OT	0331
T V _{1a}	0338

ION CHANNELS ■

Family/assay	Ref.
VOLTAGE-GATED CHANNELS	
Ca ²⁺ channels	
T L	0339
K ⁺ channels	
T K _{ATP}	0340
MEMBRANE LIGAND-GATED CHANNELS	
GABA	
T GABA _A	0314
Purinergic	
T P2X	0332
Serotonin	
T 5-HT ₃	0335

NON-KINASE ENZYMES ■

Family/assay	Ref.
PROTEASES	
Metalloproteases	
T ACE	0302
T ECE-1	0313

 tissue assay  human

■ TESTING CONDITIONS

■ SUGGESTED TESTING

Qualitative screening

- Assessment of both agonism and antagonism at 3 concentrations in duplicate (2 tissues)
- Suggested test concentrations: 1, 3 and 10 times the IC_{50}/K_i binding value

Catalog prices correspond to this combination.

Follow-up testing for quantitative determination of pharmacological parameters

- EC_{50}/pD_2 values of agonists at 6-8 concentrations in duplicate
- IC_{50} values of antagonists at 6-8 concentrations in duplicate
- pA_2 or pD'_2 values of antagonists at 3 concentrations, each at least in triplicate

Prices upon request.

■ SAMPLE SIZE (including IC_{50} follow-up studies)

For screening at 3 concentrations in duplicate (2 tissues) or follow-up: sufficient amount to prepare 150 μ l of a 1000-fold concentrated solution relative to the highest test concentration for each assay.

Maximum tolerable final DMSO concentration: 0.1%.

■ PROTOCOL

► Agonism/antagonism assessment

The agonist activity is evaluated by exposing the tissues to increasing concentrations of the compounds. Where an agonist-like response is obtained, the reference antagonist is tested against the highest test concentration to confirm the involvement of the receptor studied in this response. A reference full agonist is used as a positive control.

The antagonist activity is evaluated by testing increasing concentrations of the compounds against a single concentration of the reference agonist. A reference antagonist is used as a positive control.

► EC_{50}/pD_2 values for agonists, IC_{50} values for antagonists and maximum responses (E_{max})

Same protocols as above, with full concentration response curves.

► pA_2 or pD'_2 values for antagonists

Concentration-response curves to the reference full agonist are obtained in the absence (control) and in the presence of a fixed concentration of the compound. The resulting changes in the agonist curves are used to determine a pA_2 value for competitive antagonism (decrease in the pD_2 value of the agonist) or a pD'_2 value for non-competitive antagonism (decrease in the maximum response to the agonist).

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

■ DELIVERABLES

Mean percent of control values, EC_{50}/pD_2 , IC_{50} , pA_2 or pD'_2 values.

Concentration-response curves.

■ DATA TURNAROUND

Complete data set is typically available 4 weeks after receipt of the compound at the testing site.

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 CATALOG ONLINE](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



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