

## MULTISCREEN™ 50-GPCR+ Safety/Liability Functional Assay Panel

Multispan's panel has been specifically designed to assist researchers streamline the process of in vitro safety pharmacology profiling. With Multispan's Safety Functional Assay Panel, compounds can be rapidly evaluated while accurately evaluating potential risks from off-target effects. Waste less time on investigating suboptimal leads, enabling you to select the best drug candidates to bring to market as quickly as possible, with the MULTISCREEN™ 50-GPCR Safety Functional Assay Panel.

### Purpose

- To identify compound liabilities.
- Profiling hits, leads, and new chemotypes in a well-defined 50-GPCR assay panel implicated in central nervous system, cardiac, pulmonary, and gastrointestinal safety/liability concerns.
- Using cellular functional assays avoiding binding assay limitations for allosteric or non-functional binders and to remove binding false positives.

### Scope

- Customer may submit compounds with instructions on handling, solubility, and storage.
- Multispan shall profile compounds by FLIPR in Ca<sup>++</sup> OR cAMP assays in **agonist and antagonist modes**, single point duplicates or 10-point duplicate dose-response curves according to the customer's requirements.

### Timing

- Data will be delivered 2 weeks after receipt of compounds, P.O., and 50% upfront payment.

### Data Quality

- One control agonist shall be run for each target.
- CV < 10% and EC50 value of the control compound DRC from run to run shall not exceed 10<sup>1/2</sup>.
- Any questionable data shall be repeated at no additional charge.

Targets									
Family	Receptor	Catalog	Assay Format		Family	Receptor	Catalog	Assay Format	
			Ca <sup>++</sup>	cAMP				Ca <sup>++</sup>	cAMP
Adenosine	A1	C1427B	✓	✓	Motilin	Motilin	H1297	✓	
	A2A	C1428	✓	✓	Muscarinic	M1	C1022-1a	✓	
	A3	C1430	✓	✓		M2	C1023-1	✓	✓
Adrenergic	alpha1A	C1431-1a	✓			M3	C1024-1	✓	
	alpha1B	C1432-1	✓		M4	CG1025-1	✓	✓	
	alpha1D	C1433-1	✓		Neuropeptide Y	Y1	CG1273-1	✓	✓
	alpha2A	C1434-1	✓	✓	Opioid	delta	CG1351-1	✓	✓
	alpha2B	C1435-1	✓	✓		kappa	CG1352-1	✓	✓
	alpha2C	C1436-1	✓	✓		mu	CG1350-1	✓	✓
	beta1	beta1	C1437-1	✓	✓	Prostanoid	EP2	CG1202	✓
beta2		C1438-1a	✓	✓	EP3		CG1203-1	✓	✓
Angiotensin	AT1	H1417	✓		Purinergic	P2Y1	C1160-3	✓	
Bradykinin	B1	C1198a	✓		Serotonin	5-HT1A	C1319	✓	✓
	B2	H1199	✓	✓		5-HT1B	C1320	✓	✓
Calcitonin	CGRP	CG1515-1	✓	✓		5-HT2A	C1324-1	✓	
Cannabinoid	CB1	C1229	✓	✓		5-HT2B	C1325	✓	
Cholecystokinin	CCK1	C1038	✓			5-HT2C	C1326-1	✓	
	CCK2	H1039	✓			5-HT4	C1330	✓	✓
Dopamine	D1	H1335	✓	✓		5-HT6	C1331-1	✓	✓
	D2	C1336	✓	✓	5-HT7	C1334	✓	✓	
Histamine	H1	C1027	✓		Somatostatin	SST1	CG1345-1	✓	✓
	H2	C1028	✓	✓	Vasoactive Intestinal Peptide	VPAC1	C1292-1	✓	✓
	H3	CG1029	✓	✓		VPAC2	H1293	✓	✓
Endothelin	ETA	C1216-1	✓	✓	Vasopressin	V1A	C1042-1	✓	
	ETB	C1217-1	✓			V1B	C1043-1	✓	

\*Note: Endogenous assays are highlighted in green