



**Human Cannabinoid Type 1 Receptor  
(CB1R; CNR1)  
Reporter Assay System**

**384-well Format Assays**  
Product # IB19002

▪

**Technical Manual**  
*(version 8.0)*

**[www.indigobiosciences.com](http://www.indigobiosciences.com)**

3006 Research Drive, Suite A1, State College, PA 16801, USA

Customer Service:  
814-234-1919; FAX 814-272-0152  
[customerserv@indigobiosciences.com](mailto:customerserv@indigobiosciences.com)

Technical Service:  
814-234-1919  
[techserv@indigobiosciences.com](mailto:techserv@indigobiosciences.com)