

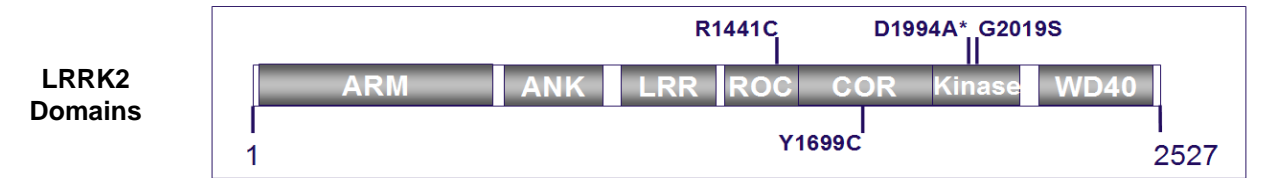
Tools for Profiling Inhibitors for Full-Length LRRK2

Justin Wetter¹, Kara Machledit¹, Ryan Fisher¹, Robert Horton¹ and David Piper², Thermo Fisher Scientific, Madison(1) and Carlsbad(2), USA

ABSTRACT

Parkinson's Disease (PD) is a progressive neurological disorder that results from degeneration of neurons in a region of the brain that controls movement, impacting an estimated 6.3 million people worldwide. To date, mutations in the leucine-rich repeat kinase 2 (LRRK2) gene are the most common genetic cause of PD. LRRK2 is a protein kinase now being aggressively pursued as a target for PD therapeutics. However, studies of LRRK2 function and inhibition have been hampered by a lack of tools to express LRRK2 in diverse cell types as well as high-throughput biochemical and cell based assays. To address these needs, we have expressed and purified a panel of full-length, disease relevant human LRRK2 isoforms, as well as BacMam particles encoding both untagged and green fluorescent protein (GFP) tagged full-length LRRK2 (WT and disease relevant mutants). This poster describes biochemical and cellular assay formats to exploit these new reagents. Together these technologies will advance the field of PD research and therapeutics by addressing the currently unmet need for modular over expression, inhibitor binding assays and cellular kinase assays for full-length LRRK2.

INTRODUCTION

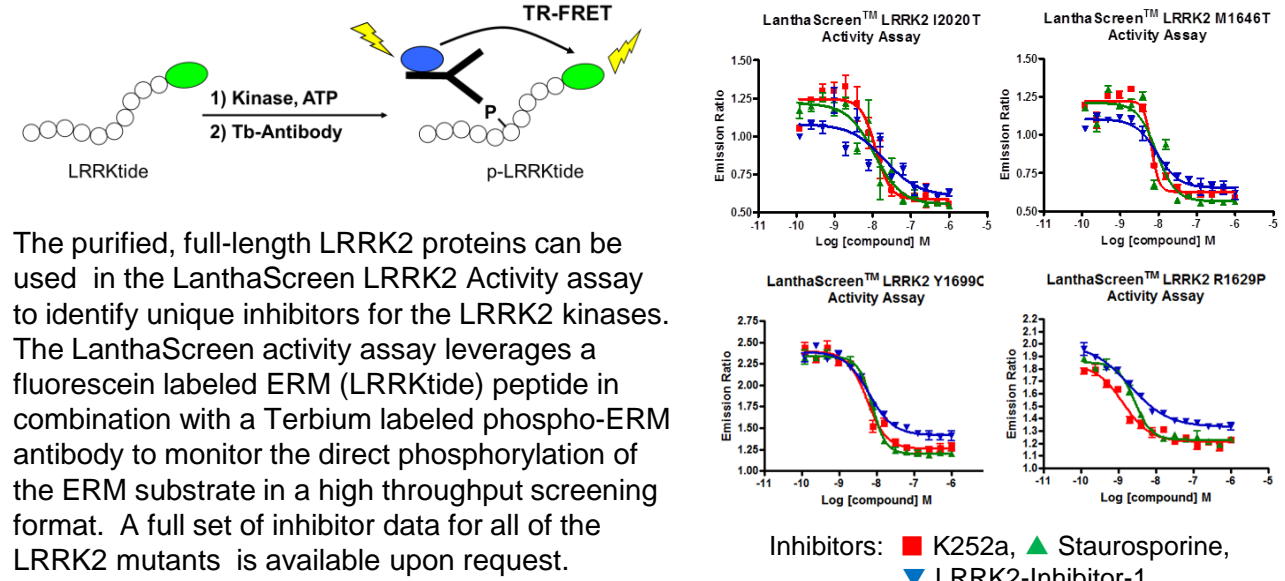


The normal function of LRRK2 is unknown but it is expressed in many tissues and organs including the brain. LRRK2 is reported to be associated with microtubules, cytoskeleton rearrangements, synaptic vesicle trafficking, neurite outgrowth and translational control. LRRK2 is a large protein (2527 amino acid) belonging to the ROCO family of proteins defined by the GTP-binding Ras of complex protein (ROC) and carboxy-terminal of Roc (COR) domains that confer weak GTPase activity. A serine/threonine protein kinase domain is positioned next to the ROC and COR domains completing the catalytic core of LRRK2. The catalytic core is flanked by leucine-rich repeats (LRR) on one side and a WD40 domain on the opposite side, both domains are thought to function as protein-protein interaction domains. Over 40 mutations of LRRK2 have been reported but only mutations in the catalytic core of LRRK2 segregate with PD in a Mendelian fashion suggesting the enzymatic functions of LRRK2 play a key role in disease pathogenesis. Mutations in the ROC GTPase domain (R1441C, R1441G, R1441H) and the COR domain (Y1699C) are reported to reduce GTPase activity in in-vitro studies. The most frequent PD associated LRRK2 mutation G2019S is located in the activation loop of the kinase domain resulting in increased kinase activity and reported to cause neurotoxicity in cultured cells.

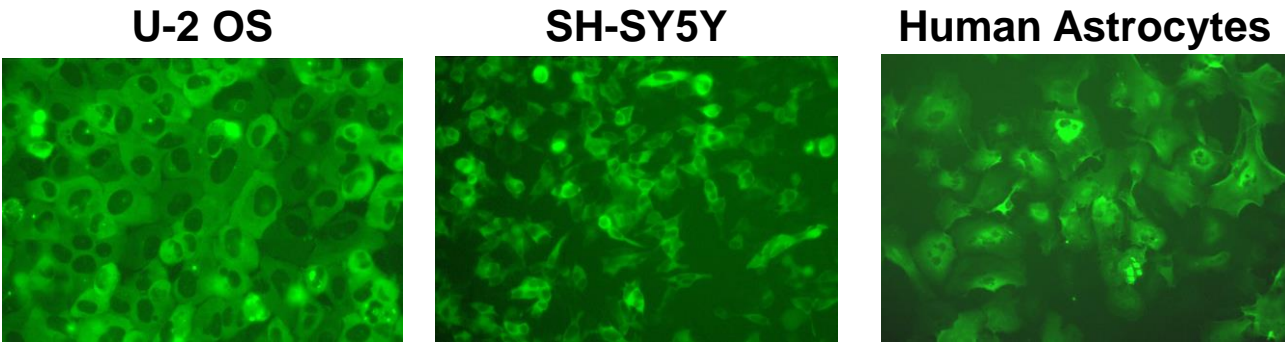
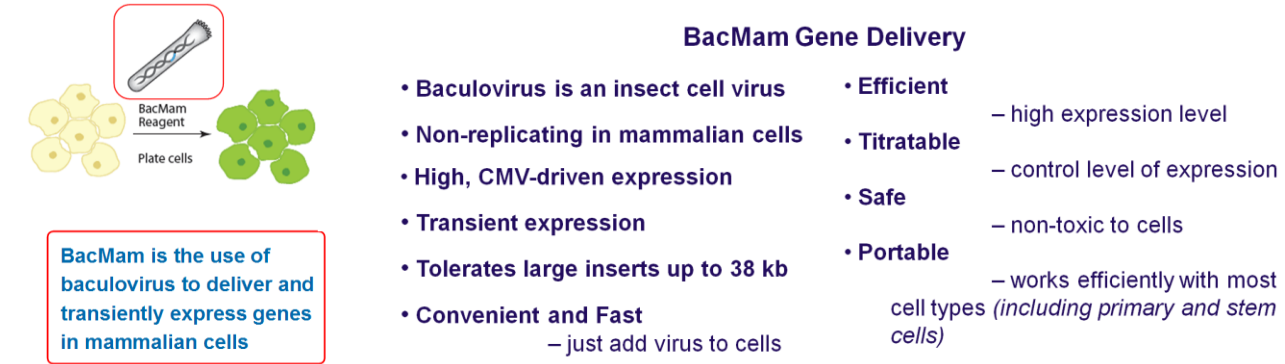
Expression and Purification of Full-Length LRRK2 protein

A BacMam expression system was successfully leveraged to express both full-length wildtype and mutant forms of LRRK2 protein in mammalian cells. The protein was purified using a DYKDDDDK-tag and tested for activity using the ERM (LRRKtide) peptide substrate (RLGRDKYKTLRQIRQ) in a radiometric assay.

LanthaScreen™ Biochemical Activity Assay for LRRK2

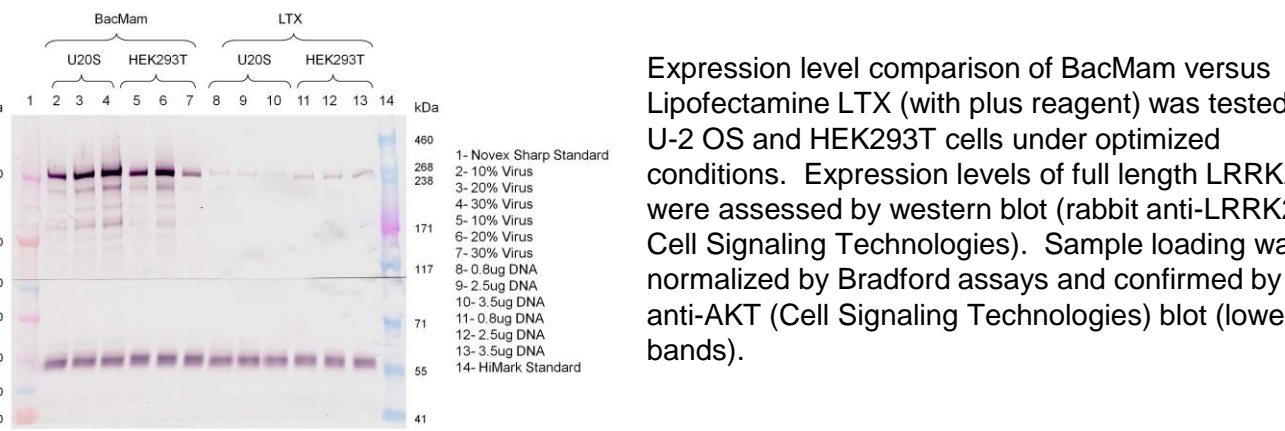


BacMam Enabled Expression of Parkinson's Disease Drug Target LRRK2

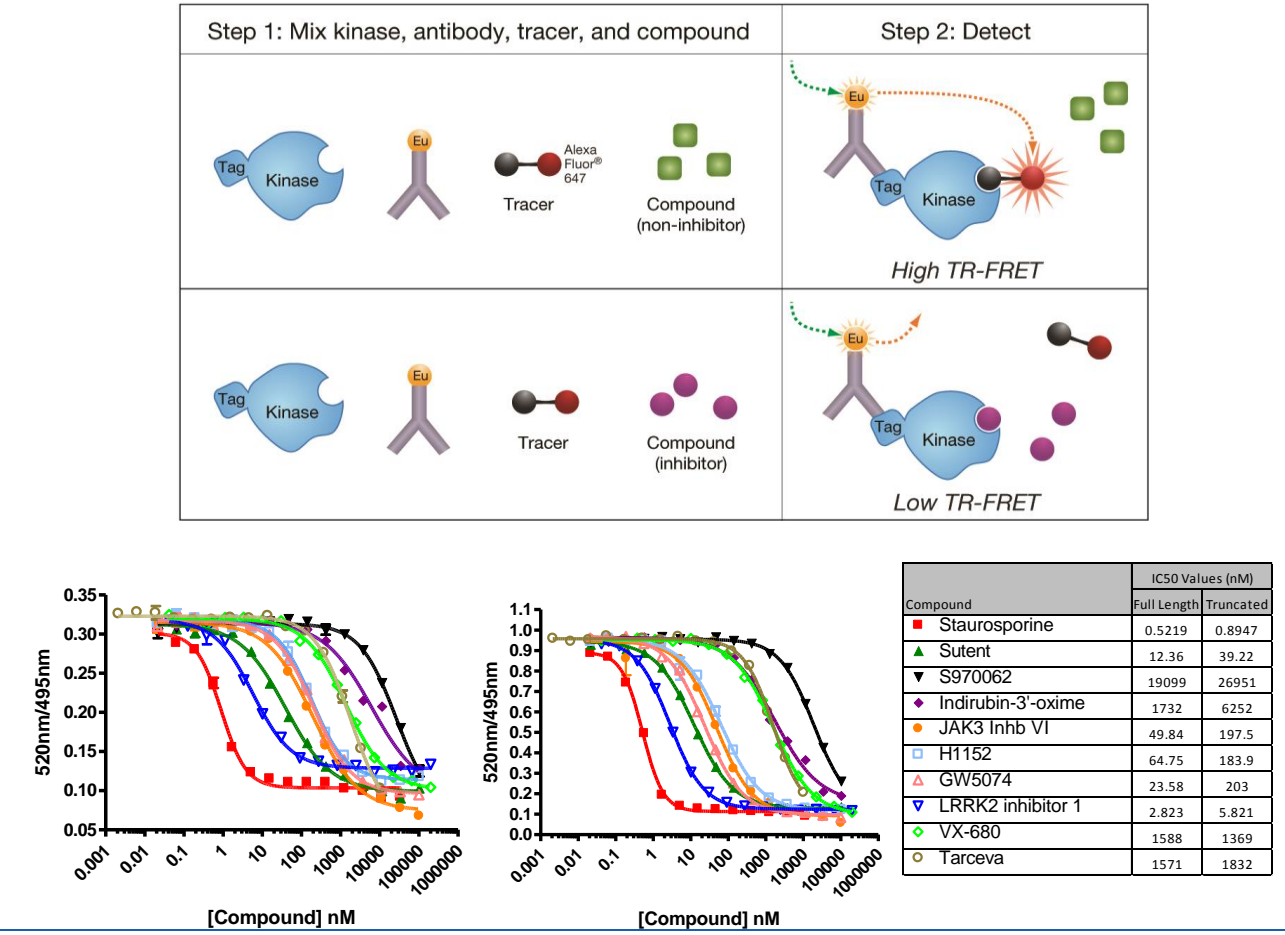


U-2 OS cells, SH-SY5Y Cells and human astrocytes (N7805100) were transduced with BacMam viruses of GFP tagged LRRK2 (A14170) and mutants. Cells were exposed to 20% virus for 24 hours, media was exchanged and cells were allowed to incubate an additional 24 hours before imaging by fluorescent microscopy

Expression Level Comparison of BacMam vs Lipid Transfection

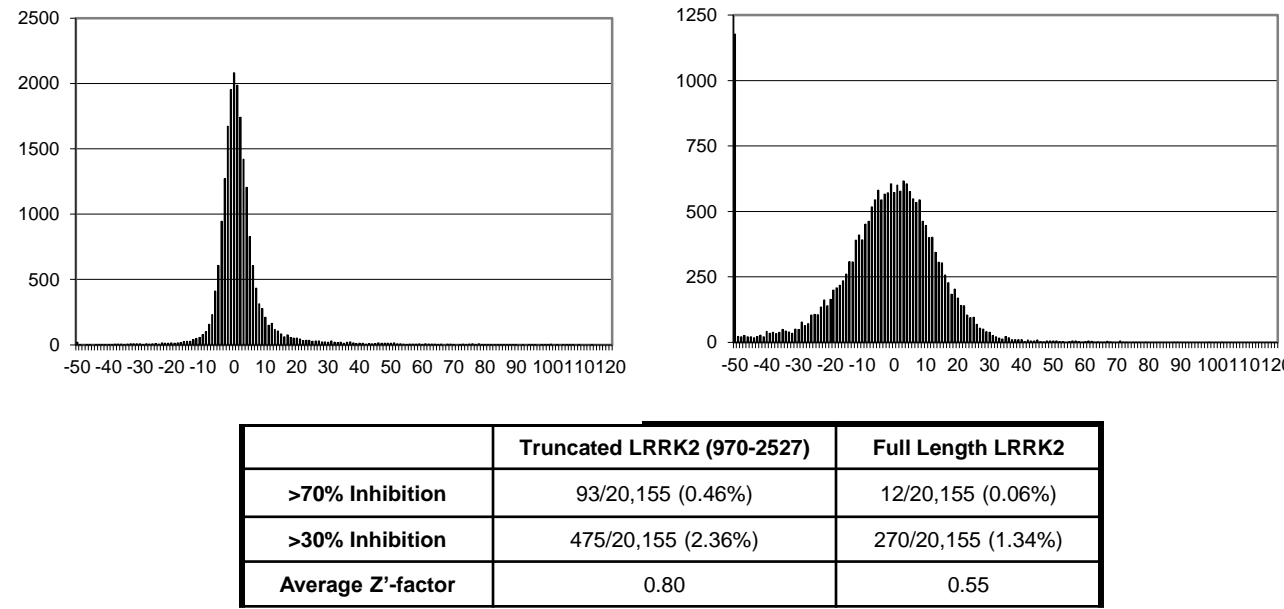


Kinase Binding Assay Utilizing Lysates from BacMam LRRK2 Transduced HEK293T Cells



BacMam LRRK2-GFP reagent was used to transduce HEK293T and generate a cellular lysate containing full length LRRK2. A lysate based Kinase Binding assay was developed using the full-length LRRK2 lysate and compared to purified truncated LRRK2. Optimal lysate and tracer concentration were determined. 5nM truncated kinase (PV4873) or 3nM full length kinase (A14171), 2 nM Eu-anti GFP antibody (A14173) and 20 nM Tracer 236 (PV5592) were used to assay binding of several kinase inhibitors to LRRK2.

Enamine Library Screen Using Full Length LRRK2 Kinase Binding Assay



Top Library Hits For LRRK2 Followed-Up Using Multiple Assay Formats

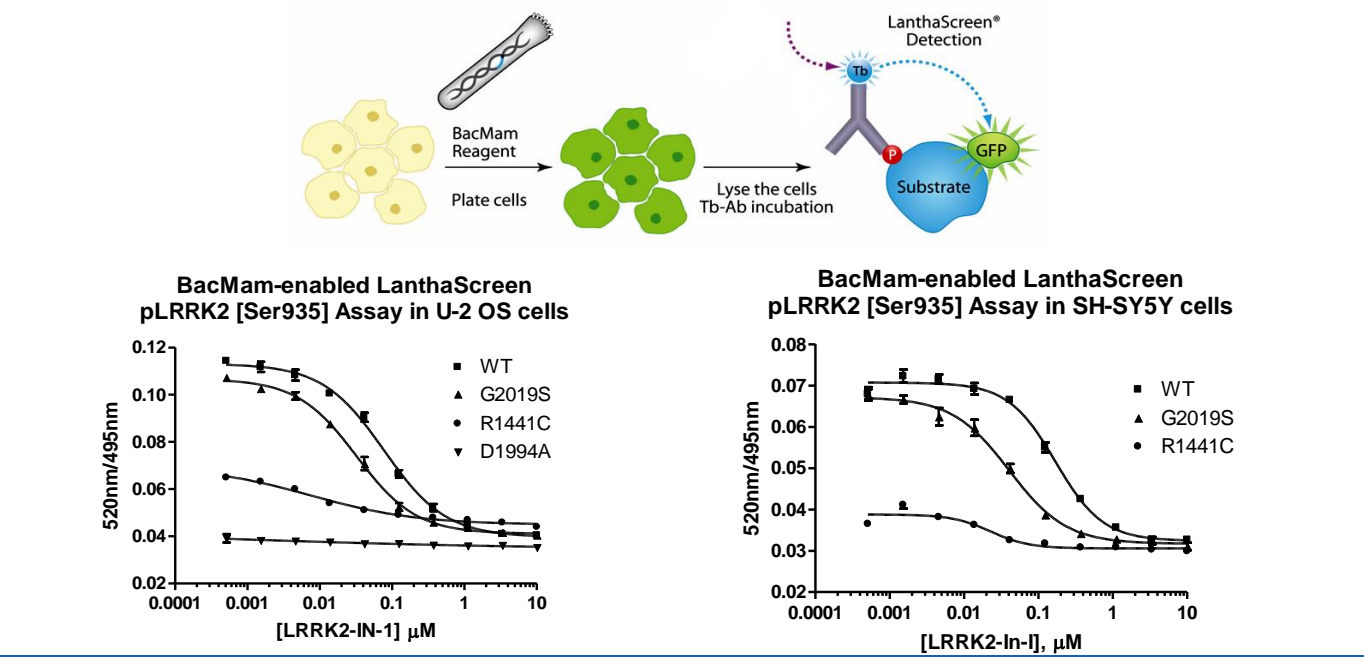
IC ₅₀ (μM)	Cellular				Biochemical Binding										Biochemical Activity	
	SH-SY5Y cells	U-2 OS cells	WT	WT	FL Lysate-based binding assay	G2019S	WT	R1441C	G2019S	WT	R1441C	D1994A	I2020T	G2019S	WT	WT
LRRK2-IN-1	0.094	0.22	0.055	0.11	0.005	0.006	0.005	0.003	0.002	0.003	0.002	0.004	0.02	0.02	nd	nd
JAK3 Inh. VI	0.73	1.2	0.21	1.4	0.23	0.2	nd	0.04	0.05	nd	nd	nd	nd	nd	nd	nd
Indirubin-3-oxime	3.8*	>10	3.9*	1.9*	9.3	6.2	nd	2.1	1.7	nd	nd	nd	nd	nd	nd	nd
Compd 2	>20	>20	>20	>20	0.13	0.15	1.9	3*	3*	5.4*	>20	>20	>20	>20	>20	>20
Compd 5	>20	>20	>20	>20	3.3	1.8	7.2	1.2	1	1.2	1.1	1.3	0.49	>20	>20	>20
Compd 6	0.53	0.76	0.2	0.31	0.09	0.06	0.06	0.025	0.012	0.017	0.011	0.029	0.18	0.18	0.18	0.18
Compd 10	>10	>10	0.37*	0.59*	2	1.9	4.2	0.09	0.11	0.21	0.11	0.23	1.3*	1.1*	1.1*	1.1*
Compd 17	>20	>20	20.9*	>20	0.84	0.8	0.59	0.39	0.26	0.34	0.24	0.52	1.8	2	2	2
Compd 25	0.17	0.88	0.99	1.22	0.03	0.18	0.12	0.01	0.019	0.028	0.014	0.053	0.11	0.26	0.26	0.26
Compd 27	nd	nd	0.44	1.29	0.25	0.274	0.6	0.054	0.034	0.1	0.056	0.19	0.54	0.88	0.88	0.88
Compd 30	>20	>20	>20	>20	8.5	1.6	1.3	10.6*	18.3*	>20	>20	11.1*	>20	>20	>20	>20

Top Library Hits Profiled Against Kinase Panel To Determine Compound Specificity

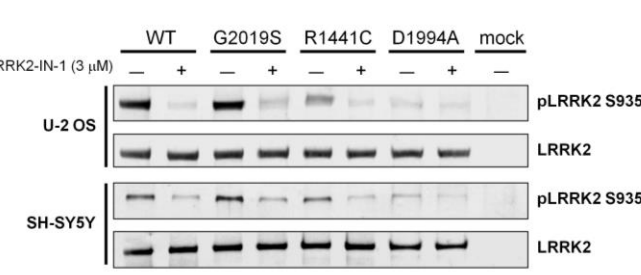
	LRRK2 IN-1	GW5074	Compd 2	Compd 5	Compd 6	Compd 10	Compd 17	Compd 25	Compd 27	Compd 30
>80% inhibition	7	126	0	216	52	8	115	14	221	5
20-80% inhibition	58	93	8	5	170	142	103	154	1	66

The lysate binding assay performed well in a high throughput screening (HTS) format when compared against the biochemical binding assay in a screen of the Enamine library - a drug like set of 20,155 compounds. Top hits from both screens were followed-up with IC₅₀ curves in multiple assay formats including lysate and biochemical binding assay, cellular assay and biochemical activity assay. Selected compounds were also profiled against a panel of 223 kinases to assess compound specificity.

Cellular Assay for Monitoring the Phosphorylation of LRRK2 Ser935



Assay Results Confirmed by Western Blot

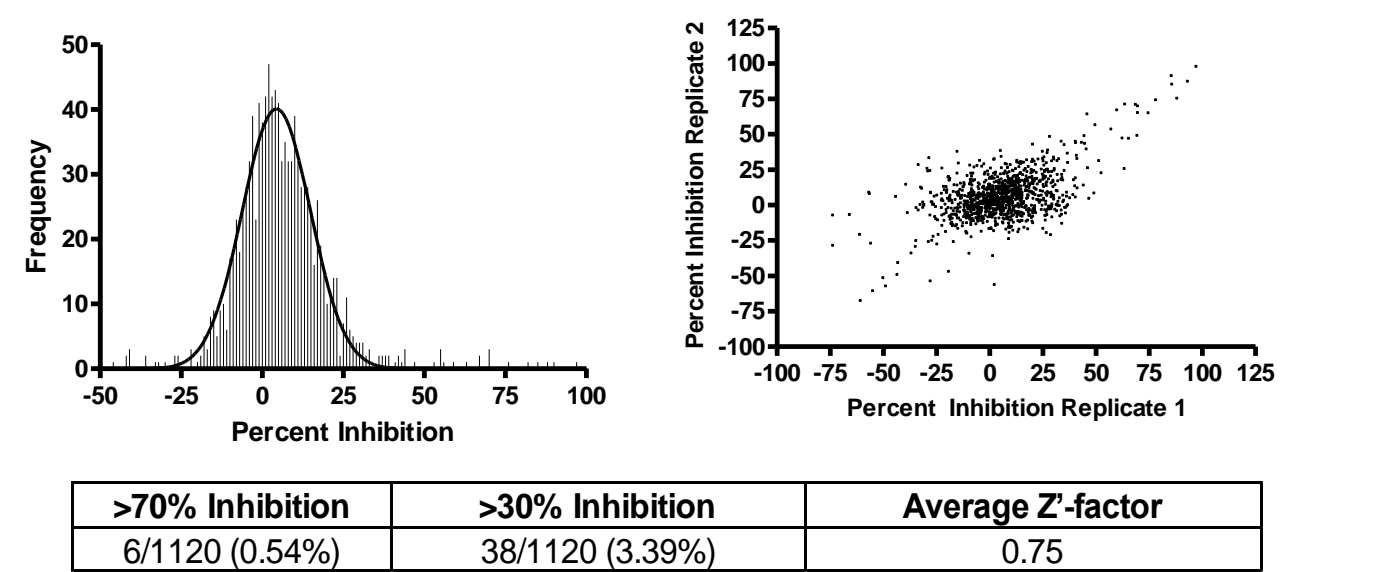


IC50 Value Summary Table

IC50 (μM)	U-2 OS		SHSY5Y		HEK293T	
Compound Name	WT	G2019S	WT	G2019S	WT	G2019S
LRRK2-IN-1	0.08	0.06	0.22	0.094	0.03	0.01
JAK3 Inh. VI	0.24	0.18	1.2	0.73	1.17	0.31
Indirubin-3-oxime	5.12	2.54	>10	3.8*	5.96	3.55
Sumitribin	0.27	0.15	0.23	0.11	0.60	0.48
GW5074	>20	0.03*	>20	>20	>20	0.28*
H-1152	2.80	0.85	1.7	0.62	3.88	1.43
H-89	>20	>20	>20	>20	nd	nd
Stauroporin	0.001	0.001	nd	nd	0.001	0.002

Using Time-Resolved Förster Resonance Energy Transfer (TR-FRET) technology, we developed a high-throughput compatible homogenous cellular assay for monitoring the LRRK2 phosphorylation at Ser935. LRRK2-GFP fusion proteins were transiently expressed in a variety of cell backgrounds via BacMam gene delivery system. Cells were plated in 384-well assay plates. The phosphorylation at Ser935 in these cells was detected using a terbium labeled anti-Ser935 phosphorylation specific antibody* that generates TR-FRET signals between terbium and GFP. Consistent with previous reports and our western blot results, wild type and G2019S LRRK2 are constitutively phosphorylated at Ser935 in cells measured by TR-FRET. The phosphorylation level was reduced for R1441C mutant* and little could be detected for the kinase activity-dead mutant D1994A*. The TR-FRET cellular assay was further validated with reported LRRK2 inhibitors including LRRK2-IN-1 and further confirmed that inhibition of LRRK2 kinase activity can reduce the phosphorylation level at Ser935.

Tocris Mini Library Screen Using LRRK2 pSer935 Cellular Assay



>70% Inhibition	>30% Inhibition	Average Z'-factor
6/1120 (0.54%)	38/1120 (3.39%)	0.75

To demonstrate the utility of the LRRK2 cellular assay in HTS applications a compound screen was performed using the Tocris mini library – a collection of 1120 biologically active compounds.

CONCLUSIONS

An extensive set of tools have been developed to enable the discovery of kinase inhibitors for PD target LRRK2. These tools allow for HTS compatible screens in multiple formats that provide detailed mechanistic data on how compounds are binding LRRK2 and inhibiting kinase activity of LRRK2 and its mutants in biochemical and cellular environments.

ACKNOWLEDGEMENTS

We would like to thank Spencer B. Hermanson, Connie Lebakken, Laurie Reichling, Ben Buscher,, and Steve Riddle for their contribution to this poster.

TRADEMARKS/LICENSING

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. For Research Use Only. Not for use in diagnostic procedures.

