

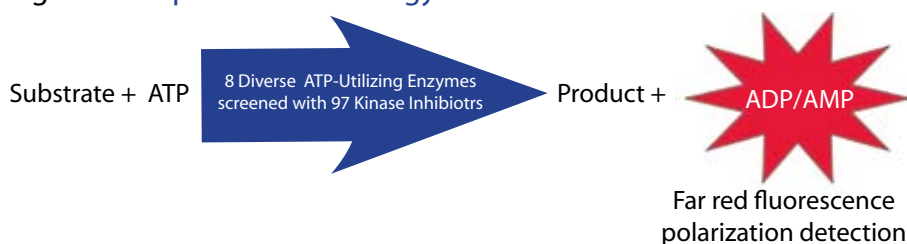
Beyond Kinases: Interrogating the Purinome

Tony A. Klink, Karen Kleman-Leyer, Matt Staeben, Robert G. Lowery
BellBrook Labs, Madison, Wisconsin, USA 53711

Introduction

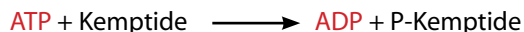
The development of ATP-site ligands as protein kinase inhibitors raises questions about the scope of their off-target effects as well as possibilities for the therapeutic targeting of other ATP-utilizing enzymes. Adenine nucleotides are interconverted by diverse proteins including, in addition to kinases, other types of transferases, phosphodiesterases, membrane transporters, DNA modifying enzymes and molecular chaperonins. Despite high diversity in the sequence motifs and folds that bind adenine nucleotides, there are commonalities in the ATP interaction networks across functionally diverse enzymes. To explore the ligand selectivity of ATP-binding sites, we screened a kinase-focused library across diverse ATP-utilizing enzymes using adenine nucleotide detection as a generic assay method. The assays rely on highly selective antibodies that distinguish between nucleotides on the basis of a single phosphate group. Homogenous fluorescent polarization assays have been developed for both ADP and AMP, making it possible to interrogate a diverse panel of otherwise intractable ATP-utilizing enzymes in an HTS format. Of the eight ATP-utilizing enzymes used in the study, we found micromolar interactions of protein kinase inhibitors with one mammalian target and one bacterial target, indicating the potential for off-target effects. The approach used provides a framework for more systematic efforts to map the ligand selectivity of ATP-utilizing enzymes. It also raises the possibility of leveraging the large body of kinase inhibitor chemoselectivity data to address other target families in the purinome.

Figure 1 Experimental Strategy

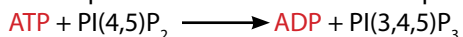


ADP-Producing Enzymes

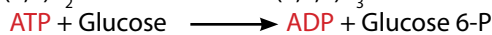
Protein Kinase A (Human):



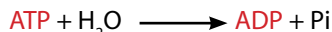
PI3 α Kinase (Human):



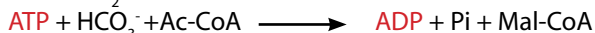
Hexokinase (Yeast):



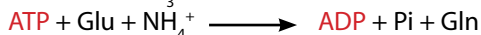
HSP73 (Bovine):



Acetyl CoA Carboxylase (Rat):



Glutamine Synthetase (M. tuberculosis):



AMP-Producing Enzymes

Ubiquitin Ligase E1 (Human):



Acyl CoA Synthetase (Pseudo. sp):

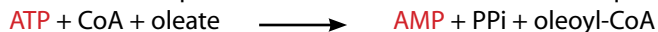
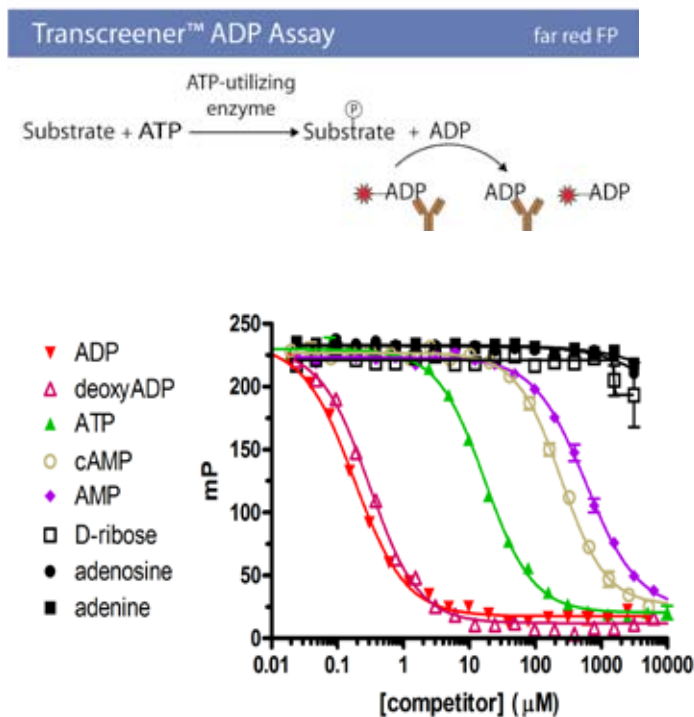
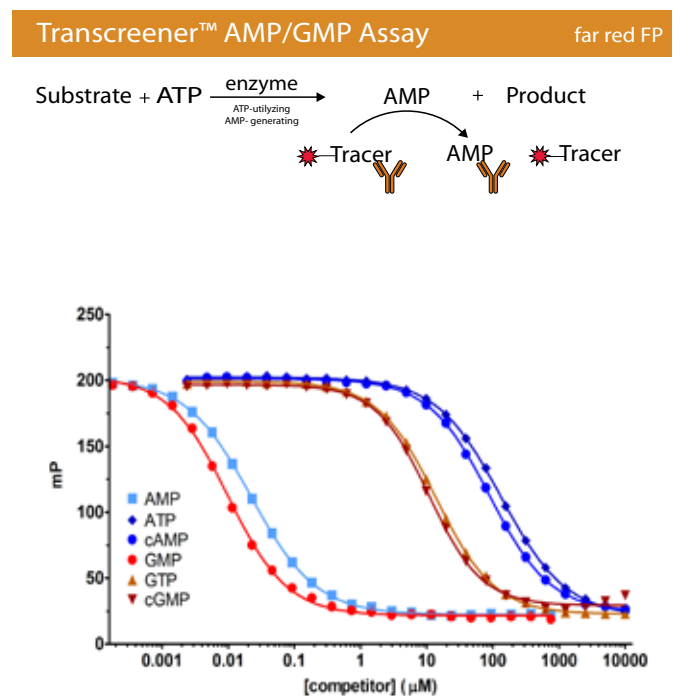


Figure 2 Transcreener Assay: Homogenous Nucleotide Detection

Transcreener™ ADP monoclonal Ab binds ADP selectively



Transcreener™ AMP/GMP Ab binds AMP selectively



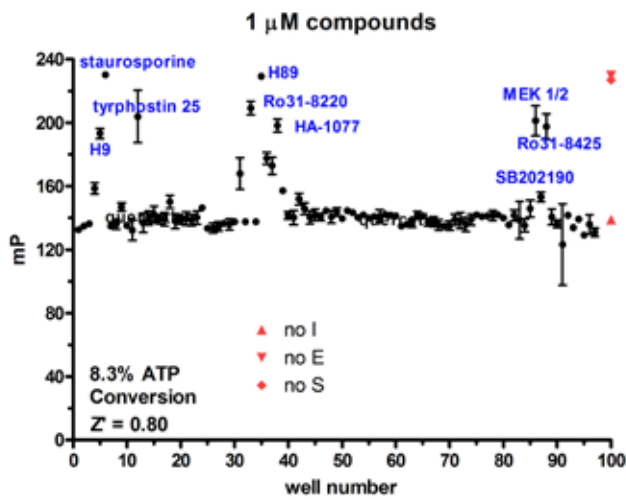
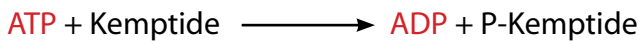
Competition curves with various adenine nucleotides demonstrate the selectivity of the Transcreener antibodies used for ADP and AMP detection. The ADP antibody discriminates ADP with greater than 100-fold selectivity versus ATP. The AMP/GMP Ab discriminates AMP with >1000-fold selectivity versus ATP or cAMP and also discriminates GMP with >1000-fold selectivity versus cGMP. The Transcreener assays are competitive fluorescence polarization immunoassays. The product of ATP-utilizing enzymes, ADP or AMP, competes with the tracer for antibody binding. Increased nucleotide concentrations result in lower polarization values.

Table 1 ATP-Utilizing Enzyme Inhibitor Profile

ATP-utilizing Enzyme	Transcreener Assay	Number of Hits (50% Inhibition at 10 μM)
PKA	ADP	19
PI3Kα	ADP	3
Acetyl CoA Carboxylase	ADP	2
Glutamine Synthetase	ADP	1
Hexokinase	ADP	0
Hsp73 ATPase	ADP	0
Acyl CoA Synthase	AMP/GMP	0
Ubiquitin E1	AMP/GMP	0

Eight ATP-Utilizing Enzymes were screened using the BIOMOL Kinase Inhibitor Library (at both 1 μM and 10 μM). In-house control compounds supplemented to increase library to 97 compounds. Micromolar inhibition seen for four of eight enzymes in the study. The number of compounds in the library that achieved 50% inhibition at 10 μM compound are listed.

Figure 3 PKA Screen



PKA reactions containing 1 μ M compound were performed in 50 mM HEPES (pH 7.5), 4 mM $MgCl_2$, 2 mM EGTA, 1% DMSO, 50 μ M Kemptide, 10 μ M ATP, and 15 ng/mL PKA (n=3). PKA control reactions (n=16) were run to 8% ATP conversion, Z' =0.8. Positive hits in the compound library screen were titrated to determine IC_{50} values.

Dose Dependency of Hits

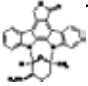
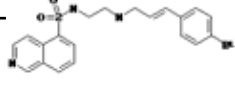
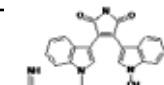
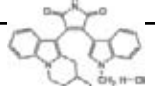
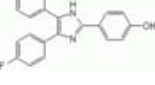
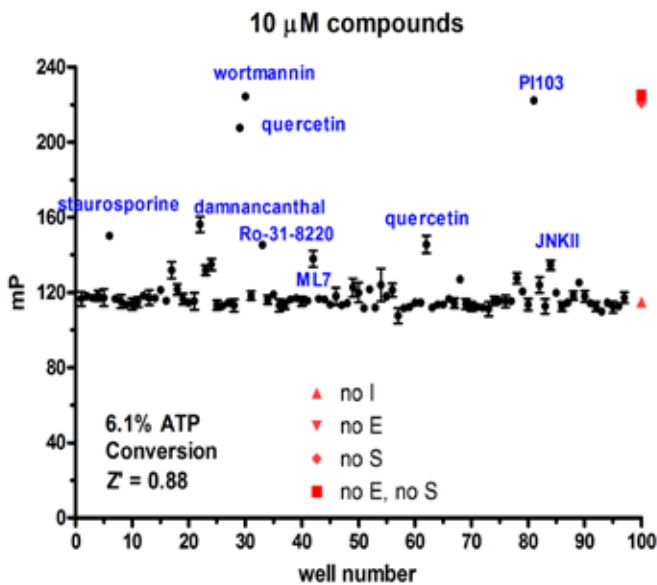
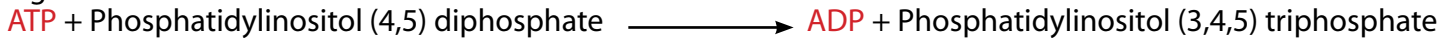
Kinase Target	Inhibitor	Structure	IC_{50}
Pan	staurosporine		4.8 nM
PKA	H89		27.3 nM
PKC/PKA	Ro-31-8220		230 nM
PKC	Ro-31-8425		451 nM
P38 MAP Kinase	SB 202190		14 μ M

Figure 4. PI3K α Screen



PI3K α reactions containing 10 μ M compound were performed in 50 mM HEPES (pH 7.1), containing NaCl (100 mM), $MgCl_2$ (4 mM), EGTA (2 mM), DTT (2 mM), DMSO (1%), ATP (30 μ M), PI(4,5) P2 C16 (30 μ M), and PI3K α (100 ng/ μ L). PI3K α control reactions (n=16) were run to 7%, Z' =0.9. Positive hits in the compound library screen were titrated to determine IC_{50} values.

Dose Dependency of Hits

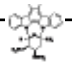
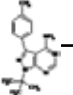
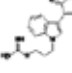
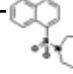
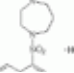
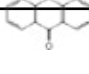
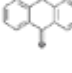
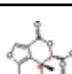
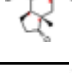

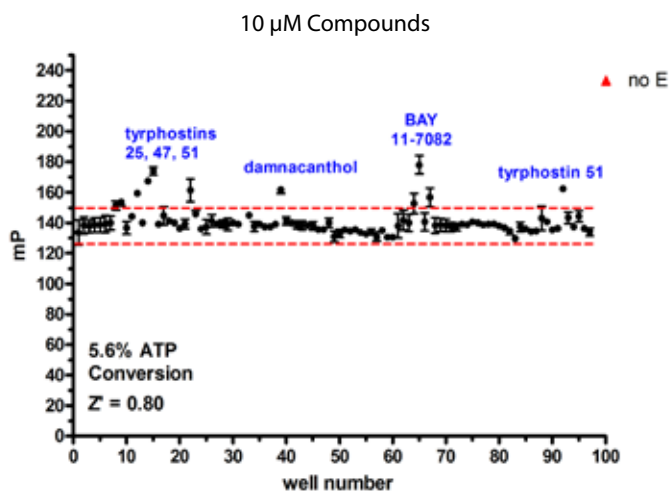
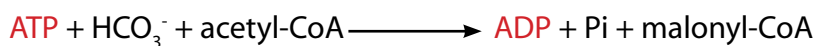
Kinase Target	Inhibitor	Structure	IC_{50}
Pan	staurosporine		9.1 μ M
Src Family	PP1		>20 μ M
PKC	Ro-31-8220		2.4 μ M
MLCK	ML7		179 μ M
MLCK	ML9		228 μ M
JNK	JNK II		>20 μ M
LCK	Damnicanthol		20 μ M
PI3K	PI 103		18 nM
PI3K	Wortmannin		14 nM
PI3K	Quercetin		>20 μ M

Figure 5. Acetyl-CoA Carboxylase Screen



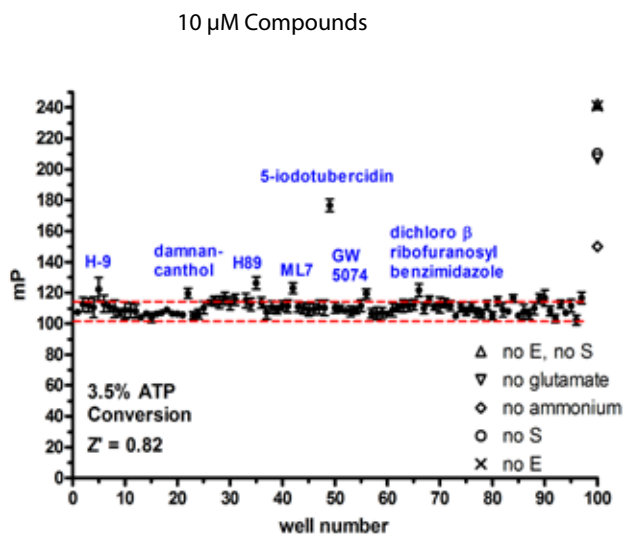
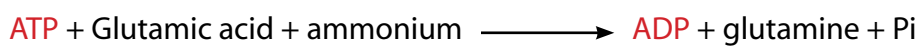
Acetyl CoA Carboxylase 1 reactions containing 10 μM compound were performed in 50 mM HEPES (pH 7.5), 4 mM MgCl_2 , 1 mM EGTA, 2 mM potassium citrate, 50 μM bicarbonate, 50 μM acetyl CoA, 1% DMSO, and 2 ng/ μL ACC1 (n=3). Acetyl CoA carboxylase control reactions (n=16) were run to 7% ATP conversion, $Z' = 0.8$.

Positive hits in the compound library screen were titrated to determine IC_{50} values.

Dose Dependency of Hits

Kinase Target	Inhibitor	Structure	IC_{50}
EGFR	Tyrphostin 25		12 μM
EGFR	Tyrphostin 47		58 μM
EGFR	Tyrphostin 51		6.8 μM
EGFR	Tyrphostin A23		9.4 μM
EGFR	Tyrphostin AG30		33 μM
EGFR	Tyrphostin B44		>100 μM
TNF α	BAY 11-7085		12.6 μM
TNF α	BAY 11-7082		6.6 μM
LCK	Damnancanthol		9.0 μM

Figure 6. Glutamine Synthetase Screen



Glutamine Synthetase (*M. tuberculosis*) reactions containing 10 μM compound were performed in 50 mM Imidazole (pH 7.1), 4 mM MgCl_2 , 0.4 mM MnCl_2 , 2 mM EGTA, 2.5 mM KCL, 1% DMSO, 0.01% Brij-35, 25 μM glutamate, 25 μM ammonium, 500 μM ATP, 800 ng/mL glutamine synthetase (n=3). Glutamine synthetase control reactions (n=16) were run to 3.5 % ATP conversion, $Z' = 0.8$. Positive hits in the compound library screen were titrated to determine IC_{50} values.

Dose Dependency of Hits

Kinase Target	Inhibitor	Structure	IC_{50}
PKA	H89		>100 μM
MLCK	ML7		>100 μM
ERK2	Iodotuber- cin		7.7 μM
cRaf	GW 5074		25 μM
CK II	Dichloro β ribofuranosyl benzimi- dazol		>100 μM
Glutamine synthetase (not in screen)	Methionine sulfoximine (not in screen)		94 μM

Materials and Methods

Materials. Glutamine Synthetase was a generous donation from Dr. Guenter Harth (UCLA) and Dr. Marcus Horwitz. (UCLA) Protein kinase A (PKA), PI3 α Kinase and ACC1 were purchased from Invitrogen (Carlsbad, CA), Millipore (Dundee, Scotland), and BlueSky (Worcester, MA), respectively. Substrates, Lipids and lipid substrates were purchased from Sigma, Avanti Polar Lipids (Alabaster, AL) or CellSignals (Columbus, OH). Inhibitors were from BIOMOL Kinase Inhibitor Library or from Sigma, Upstate, or EMD Biosciences (La Jolla, CA).

Standard plate and instrumentation settings. All assays were performed in black Corning[®] 384 Well Microplates (Corning, NY). Fluorescence intensity and polarization measurements utilizing the AMP/GMP-Alexa Fluor[®]633 or ADP-Alexa Fluor[®]633 tracers were performed on a Tecan Ultra plate reader using the following filters and settings: 612 nm excitation filter (10 nm bandwidth), 670 nm emission filter (25 nm bandwidth), 10 flashes per well, 30°C or on the Tecan Safire2[™] plate reader using the following settings: 635 nm excitation (LED), 670 nm emission (20 nm bandwidth), 10 flashes per well, 30°C.

Enzyme reactions. Enzyme concentration and reaction time were variable, but always designed to produce <10% ATP conversion. The basic assay protocol is mix and read: enzyme reactions were run in 10 μ L volumes in 384-well plates at room temperature followed by the addition of 10 μ L ADP Detection Mixture (containing EDTA and antibody/tracer complex). Reactions were equilibrated for one hour at room temperature before reading the plate.

Conclusions

- Transcreener ADP and AMP nucleotide immunodetection assays and the BIOMOL Kinase Library were used to screen eight diverse enzymes that use ATP as a substrate. The enzymes include a protein kinase, a lipid kinase, a carbohydrate kinase, a heat shock protein (ATPase), a carboxylase, a synthase, a ligase, and a microbial target involved in amino acid synthesis.
- Protein kinase inhibitors exhibit 'off target' effects for PI3K α , Acetyl-CoA Carboxylase1, and Glutamine Synthetase.
- Acetyl-CoA Carboxylase was inhibited by several tyrosine kinase inhibitors while the microbial target, Glutamine Synthetase, was inhibited by serine-threonine kinase inhibitors. PI3K α showed a mixed inhibition profile.
- The four ATP-Utilizing enzymes including Hexokinase, Hsp 73, Acyl CoA Synthase, and Ubiquitin E1 were not inhibited.

This work is supported by NIH SBIR grant CA110535-01A1.

©2008 BellBrook Labs. All rights reserved. Transcreener is a trademark of BellBrook Labs. Transcreener is a patented technology of BellBrook Labs.

BellBrook Labs 5500 Nobel Drive, Suite 250, Madison, WI 53711
866.313.7881 or 608.443.2400 www.BellBrookLabs.com