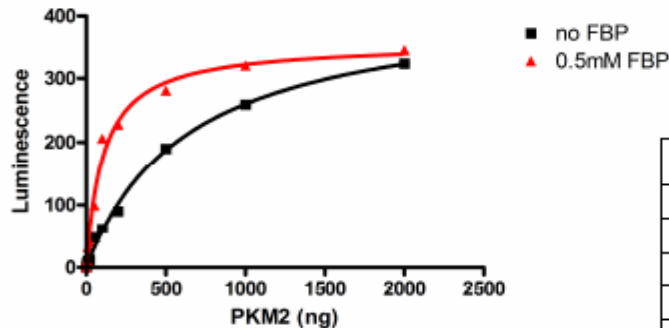


Pyruvate Kinase Profiling

Effect of FBP on PKM2 activity only at lower concentrations of PEP

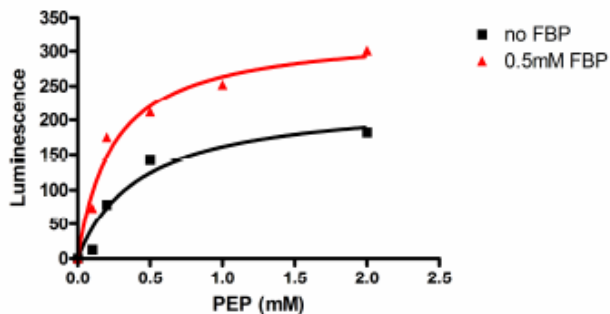


PKM2 (ng)	no FBP	0.5mM FBP
0	19	19
5	25	30
10	29	51
20	31	59
50	66	118
100	81	225
200	109	247
500	209	300
1000	278	339
2000	342	363

0.2 mM PEP and 2mM ADP in the reaction

Reference: Ikeda and Noguchi, 1998
 "Allosteric regulation of PKM2 isozyme involves a cysteine residue in the intersubunit contact"
 J Biol. Chem. Vol 273, 12227-12233, 1998.

FBP affects K_m of PEP



PEP mM	no FBP	0.5mM FBP
0	18	18
0.1	31	92
0.2	97	193
0.5	161	230
1	160	270
2	200	319

25ng PKM2 and 2mM ADP in the reaction

$K_m = 0.4224$ mM no FBP
 $K_m = 0.2467$ mM with 0.5mM FBP

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone 1.858.829.3082, Fax 1.858.481.8694
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com