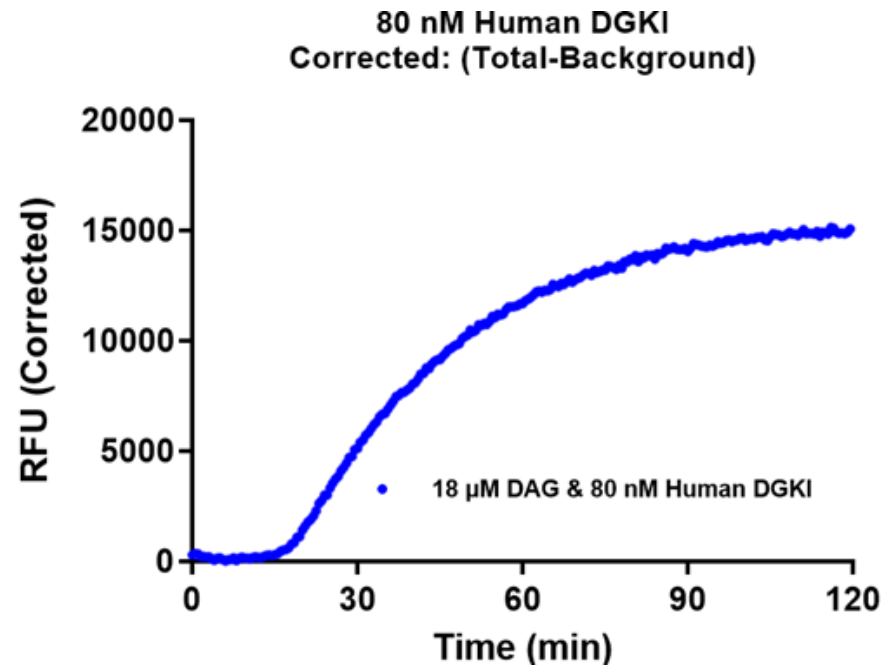
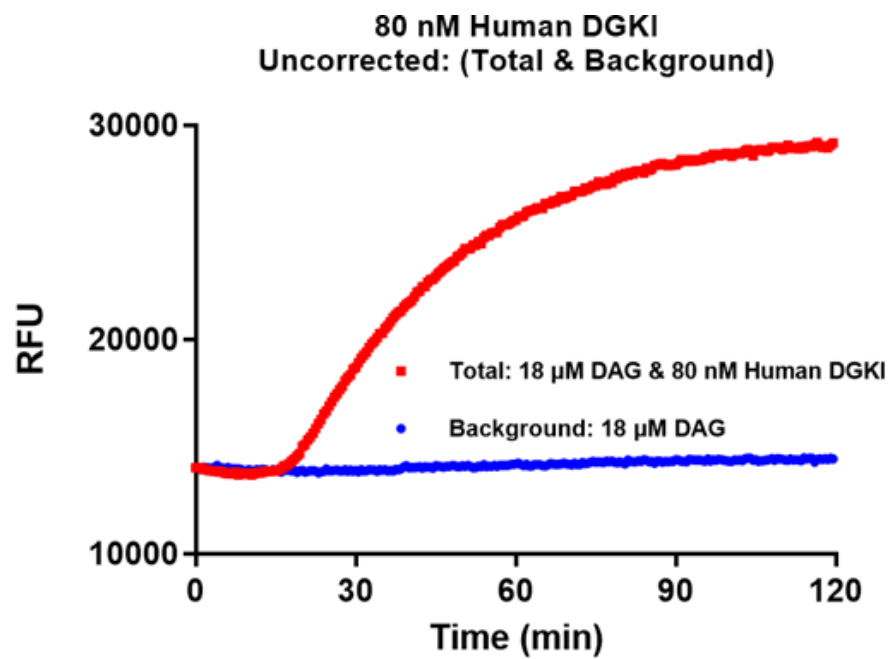


Kinase BTN-DGKI(DGKI)
Cat.# 12-407-20N
Substrate PhosphoSens[®] DAG Substrate (Assay Quant Technologies Inc.)



DAG Reaction rate: 93 ± 1 RFU/pmole/min