

PYR-41

(4-[4-(5-nitro-furan-2-ylmethylene)-3,5-dioxo-pyrazolidin-1-yl]-benzoic acid ethyl ester

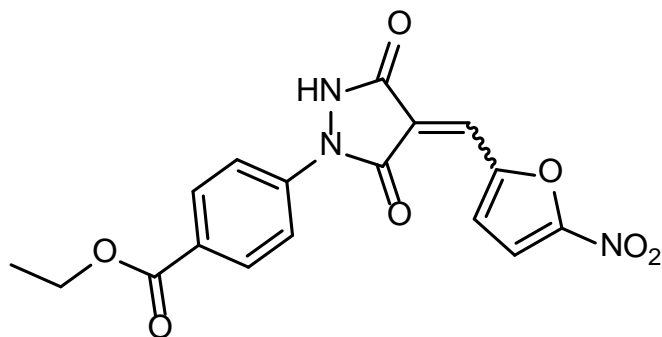
Cat. # SI9810

Background: PYR-41 is a cell permeable, irreversible inhibitor of ubiquitin activating enzyme E1 with little or no activity against E2s or E3s ($IC_{50} < 10\mu M$). It blocks degradation of p53 and inhibits NF- κ B-dependent pathways. PYR-41 also increases SUMOylation of proteins.

Application: Inhibition of ubiquitin activating enzyme E1.

Product Information

CAS No.	418805-02-4
Purity:	> 98% (TLC); NMR (Conforms)
Molecular Weight:	371.3
Physical State:	Brown powder
Quantity:	10 mg, 25 mg, 50mg
Solubility:	DMSO (25mg/ml)
Storage:	Store desiccated as supplied at 4°C for up to 2 years. Store solutions at -20°C for up to one week.



Formula: $C_{17}H_{13}N_3O_7$

References

1. Yang et al (2007) Inhibitors of ubiquitin-activating enzyme (E1), a new class of potential cancer therapeutics. *Cancer Res.* 67: 9472.
2. Brahemi et al (2010) Homology modelling of human E1 ubiquitin activating enzyme. *Lett. Drug Des. Discov.* 7: 57.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2011 LifeSensors, Inc. All Rights Reserved