

Data Sheet

N-formylkynurenine

Catalog #: 73000

Lot #: 170522

Size: 2 mg

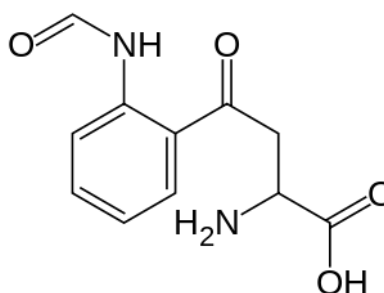
Structure:

CAS Registry #: 022-31-7

Purity: >95%

Chemical Formula: C₁₁H₁₂N₂O₄

Molecular Weight: 236.2



DESCRIPTION: N-formylkynurenine (Kyn) is an intermediate in the catabolism of tryptophan by enzymes such as indolamine 2,3-dioxygenases (IDO1 & IDO2) and tryptophan 2,3-dioxygenase (TDO). It is a formylated derivative of kynurenine and exhibits an absorption peak at 321 nm.

APPLICATIONS: Standard for measuring IDO1, IDO2, or TDO specific activity with the BPS' inhibitor screening assay kits (Cat. #72021, #72022, #72023).

STORAGE: At least 6 months when stored at -20°C as provided. *Protect from light.*

SOLUBILITY: Soluble in DMSO or methanol.

REFERENCES:

1. Seegers, N., *et al. J Biomol Screen.* 2014; **19(9)**: 1266-74.

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