

# PRODUCT DATA SHEET



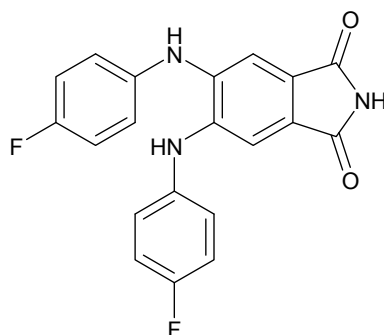
PRODUCT: DAPH-7  
CGP 53353  
4,5-bis[(4-Fluorophenyl)amino]phthalimide

CAS NO.: 145915-60-2

CATALOG NO.: EI-387

LOT NO.:

STRUCTURE:



PHYSICAL APPEARANCE: Orange solid

MOLECULAR FORMULA:  $C_{20}H_{13}F_2N_3O_2$

MOLECULAR WEIGHT: 365.3

PURITY: 98% (TLC: 50% Ethyl acetate/hexanes;  $R_f = 0.60$ )

SOLUBILITY: Soluble in: DMSO (>25 mg/mL); ethanol (5 mg/mL)

STORAGE: Store, as supplied, at  $-20^{\circ}C$  for up to 1 year. Store solutions at  $-20^{\circ}C$  for up to 3 months.

APPLICATION NOTES: DAPH-7 induces remodeling of Sup35 prion-specific intermolecular interfaces to create morphologically altered aggregates with diminished infectivity and self-templating activity. Inhibits Sup35 NM domain fibrillization with an  $IC_{50}=3.4 \mu M^1$ . DAPH-7 remodels A $\beta$ 1-42 fibers and inhibits their *de novo* assembly<sup>1</sup>. Selective PKC $\beta$ II inhibitor.  $IC_{50}$  is 410 nM for PKC $\beta$ II vs 3.8  $\mu M$  for PKC $\beta$ I.<sup>2</sup>

REFERENCES: 1. H. Wang *et al. Proc. Natl. Acad. Sci. USA* 2008 **105** 7159  
2. A. Kouroedov *et al. Circulation* 2004 **110** 91

**The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.**