

PRODUCT DATA SHEET

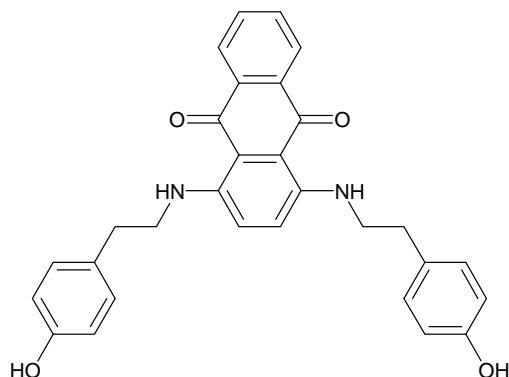


PRODUCT: BML-266
1,4-Di[(4-hydroxyphenethyl)amino]-9,10-dihydroanthracene-9,10-dione

CATALOG NO.: GR-346

LOT NO.: Q1180

STRUCTURE:



PHYSICAL APPEARANCE: Purple solid

MOLECULAR FORMULA: $C_{30}H_{26}N_2O_4$

MOLECULAR WEIGHT: 478.5

PURITY: 95% (TLC: 5% Methanol/methylene chloride; $R_f = 0.40$)

SOLUBILITY: Soluble in: DMSO (25 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: BML-266 is a structurally novel SIRT2 inhibitor ($IC_{50} = 56.7 \mu M$)¹.

REFERENCES: 1. P.H. Kiviranta *et al.* *J. Med. Chem.* 2006 **49** 7907

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.