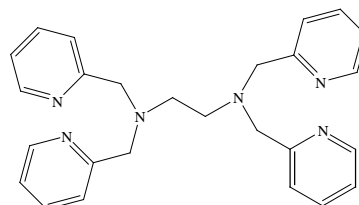


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** TPEN ( tetrakis-(2-pyridylmethyl)ethylenediamine)**CATALOG #:** 59003**MOLECULAR INFORMATION:** C<sub>26</sub>H<sub>28</sub>N<sub>6</sub>  
MW: 425**PROPERTIES:**

**Color & Form** Off-white solid  
**Purity** ≥ 95% by TLC  
**Solubility** Soluble in ethanol, DMSO or DMF

**STORAGE AND HANDLING:**

Store desiccated at ≤ 4 °C.

**APPLICATION:**

TPEN is an excellent heavy metal chelator while it does not affect Ca<sup>++</sup>, Mg<sup>++</sup>, Na<sup>+</sup> or K<sup>+</sup> concentrations. Thus, TPEN is a useful tool to identify effects of heavy metals (Zn<sup>++</sup>, Fe<sup>++</sup>/Fe<sup>+++</sup>, Cu<sup>++</sup>, and Mn<sup>++</sup>, etc.) on fluorescent indicators of Ca<sup>++</sup>, Mg<sup>++</sup>, Na<sup>+</sup>, and K<sup>+</sup>.

Ref. *J. Biol. Chem.* 260, 2719(1985).**TOXICITY:**

Unknown.

**FIRST AID:**

Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.

**Disclaimer:** Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.