

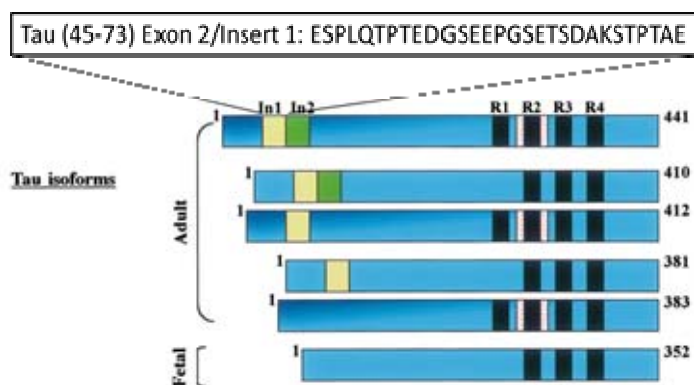


## Product Data Sheet

|                          |  |                               |
|--------------------------|--|-------------------------------|
| <b>Product Name:</b>     | Tau Peptide (45-73) (Exon 2/Insert 1 domain)   |                               |
| <b>Catalog Number:</b>   | AS-65431-1 (1 mg); AS-65431-5 (5 mg)   | Lot Number: See label on vial |
| <b>Sequence:</b>         | Glu-Ser-Pro-Leu-Gln-Thr-Pro-Thr-Glu-Asp-Gly-Ser-Glu-Glu-Pro-Gly-Ser-Glu-Thr-Ser-Asp-Ala-Lys-Ser-Thr-Pro-Thr-Ala-Glu-OH (3-letter code) |                               |
|                          | ESPLQPTEDGSEEPGSETSDAKSTPTAE (1-letter code)   |                               |
| <b>Molecular Weight:</b> | 2977.97  |                               |
| <b>Peptide Purity:</b>   | Peak Area by HPLC ≥95%   |                               |

**Storage:** Tau Peptide (45-73) (Exon 2/Insert 1 domain) is shipped at ambient temperature. Upon receipt, store lyophilized peptide at  $-20^{\circ}\text{C}$  or lower. Reconstituted peptide can be aliquoted and stored at  $-20^{\circ}\text{C}$  or lower.

**Description:** TAU proteins belong to the microtubule-associated protein (MAP) family and are involved in the pathogenesis of Alzheimer's disease. In the human brain, there are six TAU isoforms ranging from 352 to 441 amino acids in length. These isoforms vary at the carboxyl terminal according to the presence of either three repeat (3R) or four repeat (R4) domains, in addition to the presence or absence of one or two insert domains at the amino-terminus (see figure below). Tau Peptide (45-73) is a 29-amino acid long peptide derived from the Exon 2/Insert 1 domain.



### Related Products:

- Tau Peptide (74-102) (Exon 3/Insert 2 domain), Cat# AS-65432-1 (1 mg) & AS-65432-5 (5 mg)
- Tau Peptide (244-274) (Repeat 1 domain), Cat# AS-65433-1 (1 mg) & AS-65433-5 (5 mg)
- Tau Peptide (275-305) (Repeat 2 domain), Cat# AS-65434-1 (1 mg) & AS-65434-5 (5 mg)
- Tau Peptide (306-336) (Repeat 3 domain), Cat# AS-65435-1 (1 mg) & AS-65435-5 (5 mg)
- Tau Peptide (337-368) (Repeat 4 domain), Cat# AS-65436-1 (1 mg) & AS-65436-5 (5 mg)

*For Research Use Only*