



## Product Data Sheet

---

<b>Product Name:</b>	AggreSure™ beta-Amyloid (1-42), human
<b>Catalog Number:</b>	AS-72216 (0.25 mg, net peptide)
<b>Sequence:</b>	H - Asp - Ala - Glu - Phe - Arg - His - Asp - Ser - Gly - Tyr - Glu - Val - His - His - Gln - Lys - Leu - Val - Phe - Phe - Ala - Glu - Asp - Val - Gly - Ser - Asn - Lys - Gly - Ala - Ile - Ile - Gly - Leu - Met - Val - Gly - Gly - Val - Val - Ile - Ala - OH (3-letter code)  DAEFRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter code)
<b>Molecular Weight:</b>	4514.1
<b>Peptide Purity:</b>	>95%
<b>Appearance:</b>	Lyophilized white powder
<b>Peptide Reconstitution:</b>	Reconstitute peptide in either 50mM Tris/150mM NaCl (pH 7.2) or 20mM HEPES/150mM NaCl (pH 7.2) at 0.25 mg/ml. Use water bath sonicator to completely dissolve peptide if necessary. <b>Do not vortex!</b>
<b>Storage:</b>	Upon receipt, store lyophilized peptide at -20°C or lower. Do not freeze-thaw reconstituted peptide.
<b>Description:</b>	AggreSure™ beta-Amyloid (1-42) peptide was pretreated and tested for aggregation using SensoLyte® ThT Aβ42 Aggregation kit (Cat. AS-72214).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
AggreSure™ beta-Amyloid (1-40), human	72215	0.25 mg
AggreSure™ [Pyr3] β-Amyloid (3-42), human	72277	0.25 mg
Beta-Amyloid (1-40), human	24236	1 mg
Beta-Amyloid (1-42), human	20276	1 mg
[Pyr3]-beta-Amyloid (3-42)	29907	1 mg
SensoLyte® Thioflavin T Aβ40 Aggregation Kit	72213	1 kit
SensoLyte® Thioflavin T Aβ42 Aggregation Kit	72214	1 kit
SensoLyte® Thioflavin T [Pyr3] β-Amyloid (3-42) Aggregation Kit	72276	1 kit