



Product Data Sheet

Product Name:	AggreSure [Pyr3] β -Amyloid (3-42), human
Catalog Number:	AS-72277 (0.25 mg, net peptide)
Sequence:	Pyr-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) Pyr-FRHDSGYEVHHQKLVFFAEDVGSNKGAIIGLMVGGVVIA (1-letter code)
Molecular Weight:	4310.2
Peptide Purity:	>95%
Appearance:	Lyophilized white powder
Peptide Reconstitution:	To reconstitute peptide, use <i>cold</i> solvents. Add 100 μl of 1%NH₄OH, then add 900 μl of 50mM Tris/150mM NaCl (pH 7.2) to achieve a concentration of 0.25 mg/ml. Use water bath sonicator to completely dissolve peptide if necessary. <i>Do not vortex!</i>
Storage:	AggreSure [Pyr3] β -Amyloid (3-42), human is shipped at ambient temperature. Upon receipt, store lyophilized peptide at -20°C or lower. Do not freeze-thaw reconstituted peptide.
Description:	AggreSure [Pyr3] β -Amyloid (3-42), human peptide was pretreated and tested for aggregation using SensoLyte® Thioflavin T [Pyr3] β -Amyloid (3-42) Aggregation Kit (Cat. AS-72276).

Related Products:

Name	Cat #	Size
AggreSure™ beta-Amyloid (1-40), human	AS-72215	0.25 mg
AggreSure™ beta-Amyloid (1-42), human	AS-72216	0.25 mg
[Pyr3]-beta-Amyloid (3-42)	AS-29907	1mg
Beta-Amyloid (1-40), human	AS-24236	1 mg
Beta-Amyloid (1-42), human	AS-20276	1 mg
SensoLyte® Thioflavin T β -amyloid (1-40) Aggregation Kit	AS-72213	1 kit
SensoLyte® Thioflavin T β -amyloid (1-42) Aggregation Kit	AS-72214	1 kit
SensoLyte® Thioflavin T [Pyr3] β -Amyloid (3-42) Aggregation Kit	AS-72276	1 kit