

Neuroscience Proteomics Portfolio

AnaSpec, a subsidiary of Eurogentec, is dedicated to helping scientists advance neurodegenerative disease research by offering a specialized grouping of **peptides, assay kits, recombinant proteins, and dyes**. These **off-the-shelf** catalog products represent over 30 years of product innovation, and are organized by the following research areas for your convenience: **Alzheimer's disease, Parkinson's disease and Multiple sclerosis**.



Alzheimer's disease research products

Assay kits

Catalog#	Kit	Assay type	Detection
AS-72085	Sensolyte® 520 TACE (α-Secretase) Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-71144	Sensolyte® 520 β-Secretase (BACE1) Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72225	Sensolyte® 520 BACE2 Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72213	Sensolyte® Thioflavin T β-Amyloid (1-40) Aggregation Kit	Fluorimetric Assay	Fluorescent
AS-72214	Sensolyte® Thioflavin T β-Amyloid (1-42) Aggregation Kit	Fluorimetric Assay	Fluorescent
AS-72243	Sensolyte® 520 Endothelin Converting Enzyme (ECE) Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72223	Sensolyte® 520 Neprilysin Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72230	Sensolyte® Green Glutamyl Cyclase Activity Assay Kit *Fluorimetric*	Fluorogenic substrate Assay	Fluorescent
AS-72231	Sensolyte® 520 Insulin Degrading Enzyme (IDE) Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72242	Sensolyte® 520 Acetylcholinesterase Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72240	Sensolyte® Green Peptidyl Prolyl Isomerase (PINT) Activity Assay Kit *Fluorimetric*	Fluorogenic substrate Assay	Fluorescent

Dyes

Catalog#	Dye	Size
AS-88300	BSB [(trans,trans)-1-Bromo-2,5-bis-(3-hydroxycarbonyl-4-hydroxy)styryl]benzene]	5 mg
AS-83016	Congo Red *UltraPure Grade*	1 g
AS-88306	Thioflavin T *UltraPure Grade*	1 g

Recombinant Proteins

Catalog#	Recombinant Protein	Size
AS-55556-50	Tau (Tau-441), human recombinant	50 µg
AS-55556-100	Tau (Tau-441), human recombinant	100 µg
AS-55557-50	Tau (Tau-441), human recombinant, GST tagged	50 µg
AS-55557-100	Tau (Tau-441), human recombinant, GST tagged	100 µg

Peptides

β-Amyloid 1-42

Catalog#	Peptide	Size	Catalog#
AS-24224	Beta-Amyloid (1-42)	0.5mg	AS-24235
AS-20276		1mg	AS-24236
AS-20276-5		5mg	AS-24236-5
AS-20276-25		25mg	
AS-64129-05	Beta-Amyloid (1-42) • HFIP ¹	0.5mg	AS-64128-05
AS-64129-1		1mg	AS-64128-1
AS-72216	AggreSure Beta-Amyloid (1-42), human ²	0.25mg	AS-72215
AS-GMP-20276-1	GMP Beta-Amyloid (1-42), human ³	1mg	AS-GMP-24236-1
AS-21793	Beta-Amyloid (1-42) • HCl	1mg	AS-20698
AS-25383	Scrambled-beta-Amyloid (1-42)	1mg	AS-24626
AS-27276	Beta-Amyloid (42-1)	1mg	AS-22818
AS-25231	Beta-Amyloid Peptide (1-42), mouse, rat	1mg	AS-25230
AS-23524-01	Biotin-beta-Amyloid (1-42)	0.1mg	AS-23512-01
AS-24641-01	Biotin-LC-beta-Amyloid (1-42)	0.1mg	AS-24645-01
AS-23538	Cys-beta-Amyloid (1-42)	1mg	AS-23520
AS-61484-01	Beta-Amyloid (1-42)-Lys(Biotin)-NH2	0.1mg	AS-61483-01
AS-23526-01	Beta-Amyloid (1-42), FAM-labeled	0.1mg	AS-23514-01
AS-60476	Beta-Amyloid (1-42), TAMRA-labeled	0.1mg	AS-60488-01
AS-60479-01	Beta-Amyloid (1-42), HiLyte™ Fluor 488-labeled	0.1mg	AS-60491-01

β-Amyloid 1-40 equivalent

β-Amyloid 1-42

Catalog#	Peptide	Size	Catalog#
AS-60480-01	Beta-Amyloid (1-42), HiLyte™ Fluor 555-labeled	0.1mg	AS-60492-01
AS-64161	Beta-Amyloid (1-42), HiLyte™ Fluor 647-labeled	0.1mg	AS-60493
AS-63323	[Arg6]-beta-Amyloid (1-42), English Mutation	0.5 mg	AS-63321
AS-63324	[Asn7]-beta-Amyloid (1-42), Tottori-Japanese Mutation	0.5 mg	AS-63320
AS-63704	[Gly21]-beta-Amyloid (1-42), Dutch Mutation	0.5 mg	AS-62150
AS-61967-05	[Gly22]-beta-Amyloid (1-42), Arctic Mutation	0.5 mg	AS-61262
AS-62142	[Gln22]-beta-Amyloid (1-42), Dutch mutation	0.5 mg	AS-61261
AS-62148	[Lys22]-beta-Amyloid (1-42), Italian Mutation	0.5 mg	AS-62147
AS-63705	[Asn23]-beta-amyloid (1-42), Iowa Mutation	0.5 mg	AS-62145
AS-63672-1	[Cys26]-beta-Amyloid (1-42)	1 mg	AS-63671-1
AS-63738	ClearPoint™ beta-Amyloid (1-42), 13C,15N-labeled on R+K	50 µg	AS-63737
AS-29909-1	Beta-Amyloid (2-42)	1 mg	AS-29905-1
AS-63715-01	Beta-Amyloid (3-42)	0.1 mg	AS-61029
AS-29907-1	[Pyr3]-beta-Amyloid (3-42)	1 mg	AS-29906-1
AS-63317	Beta-Amyloid (11-42)	1 mg	AS-60017-1
AS-29903-1	[Pyr11]-beta-Amyloid (11-42)	1 mg	AS-29904-1

β-Amyloid 1-40 equivalent

1 HFIP treated to remove any preexisting structures for controlled aggregation studies — 2 Proprietary pre-treatment to guarantee aggregation — 3 GMP controlled production

Other Amyloid peptide fragments and related peptides

Catalog#	Peptide	Size	Catalog#	Peptide	Size	Catalog#	Peptide	Size
AS-24232	Beta-Amyloid (1-28)	1 mg	AS-60252-05	Beta-Amyloid (1-36)	0.5 mg	AS-61971	Beta-Amyloid (25-35), scrambled	1 mg
AS-62465	Beta-Amyloid (1-28)-Lys(Biotin)-NH2	0.5 mg	AS-60251-1	Beta-Amyloid (1-37)	1 mg	AS-62451	Biotin-beta-Amyloid (25-35)	1 mg
AS-24230	Beta-Amyloid (12-28)	1 mg	AS-24234	Beta-Amyloid (1-38)	1 mg	AS-62476	Beta-Amyloid (1-38), mouse, rat	1 mg
AS-65113	[Pro18]-beta-Amyloid (12-28)	1 mg	AS-65220	ClearPoint™ beta-Amyloid (1-38), 13C-Phe & Ile	0.05 mg	AS-65431-1	Tau peptide (45-73) (Exon2/Insert1 domain)	1 mg
AS-61802	Beta-Amyloid (13-28)	1 mg	AS-24296	Beta-Amyloid (1-39)	1 mg	AS-65433-1	Tau peptide (244-274) (Repeat1 domain)	1 mg
AS-62447	Beta-Amyloid (17-28)	1 mg	AS-25357	Beta-Amyloid (1-43)	1 mg	AS-65434-1	Tau peptide (275-305) (Repeat2 domain)	1 mg
AS-61970	Beta-Amyloid (1-9)	1 mg	AS-63335	Beta-Amyloid (1-55)	0.1 mg	AS-65435-1	Tau peptide (306-336) (Repeat3 domain)	1 mg
AS-64478	Beta-Amyloid (1-10)	1 mg	AS-64793	Beta-Amyloid (3-16)	1 mg	AS-65436-1	Tau peptide (337-368) (Repeat4 domain)	1 mg
AS-22819	Beta-Amyloid (1-11)	1 mg	AS-25365	Beta-Amyloid (10-20)	1 mg	AS-65409-1	Ser(OGlcNAc)400-KWK-Tau (388-411)-KK(Biotin)-NH2	1 mg
AS-62967	Beta-Amyloid (1-12)	1 mg	AS-62885	Beta-Amyloid (10-26)	1 mg	AS-60886	Humanin, HN	1 mg
AS-64480	Beta-Amyloid (1-13)	1 mg	AS-63745	Beta-Amyloid (10-35)-Lys(Biotin)-NH2	0.1 mg	AS-60887	Humanin S14G, sHNG	1 mg
AS-61798	Beta-Amyloid (1-15)	1 mg	AS-63819	Beta-Amyloid (11-17)	1 mg	AS-60888	Humanin C8A, sHNA	1 mg
AS-24225	Beta-Amyloid (1-16)	1 mg	AS-23956	Beta-Amyloid (11-22)	1 mg	AS-61930	Colivelin	1 mg
AS-62136	Beta-Amyloid (1-16)-Lys(Biotin)-LC)-NH2	1 mg	AS-62884	Beta-Amyloid (11-25)	1 mg	AS-65579	Beta-Amyloid Bri (1-23)	1 mg
AS-62002-05	Beta-Amyloid (1-15)-Lys16, HiLyte™ Fluor 488-labeled	0.5 mg	AS-62907	Biotin-LC-beta-Amyloid (15-25)	1 mg	AS-65580	Beta-Amyloid Bri (1-34)	1 mg
AS-61955	Beta-Amyloid (1-17)	1 mg	AS-62897	Beta-Amyloid (16-22)	1 mg	AS-65575	Neurogranin (48-76), human	1 mg
AS-62473	Beta-Amyloid (1-17)-Cys	1 mg	AS-61978	Beta-Amyloid (17-24)	1 mg	AS-65576	Biotin-Neurogranin (48-76), human	1 mg
AS-62966	Beta-Amyloid (1-22)	1 mg	AS-22823	Beta-Amyloid (22-35)	1 mg	AS-65577	Biotin-Neurogranin (48-76), mouse	1 mg
AS-60250-1	Beta-Amyloid (1-33)	1 mg	AS-62455	Biotin-LC-beta-Amyloid (22-41)	1 mg	AS-65578	Neurogranin (48-76), mouse	1 mg
AS-61799	Beta-Amyloid (1-34)	0.5 mg	AS-24227	Beta-Amyloid (25-35)	1 mg			
			AS-63308	Beta-Amyloid (25-35), HiLyte™ Fluor 488-labeled	0.1 mg			



Parkinson's disease research products

Recombinant Proteins

Catalog#	Recombinant Protein	Size
AS-55457	Human recombinant α-Synuclein (1-140), HiLyte™ Fluor 488 labeled	200 µg
AS-55581	Human recombinant α-Synuclein (1-140), biotin labeled	200 µg
AS-55555-100	EndoClear™ human recombinant α-Synuclein protein	100 µg
AS-55555-500	EndoClear™ human recombinant α-Synuclein protein	500 µg
AS-55555-1000	EndoClear™ human recombinant α-Synuclein protein	1000 µg
AS-56082-100	EndoClear™ mouse recombinant α-Synuclein protein	100 µg
AS-56082-500	EndoClear™ mouse recombinant α-Synuclein protein	500 µg

Catalog#	Recombinant Protein	Size
AS-56082-1000	EndoClear™ mouse recombinant α-Synuclein protein	1000 µg
AS-56083-100	EndoClear™ rat recombinant α-Synuclein protein	100 µg
AS-56083-500	EndoClear™ rat recombinant α-Synuclein protein	500 µg
AS-56083-1000	EndoClear™ rat recombinant α-Synuclein protein	1000 µg
AS-56081-100	EndoClear™ Plus human recombinant α-Synuclein protein	100 µg
AS-56081-500	EndoClear™ Plus human recombinant α-Synuclein protein	500 µg
AS-56086-100	EndoClear™ Plus mouse recombinant α-Synuclein protein	100 µg
AS-56086-500	EndoClear™ Plus mouse recombinant α-Synuclein protein	500 µg



Multiple Sclerosis research products

Assay kits

Catalog#	Kit	Assay type	Detection
AS-54465	Sensolyte® Anti-MOG(35-55) IgG Quantitative ELISA Kit (mouse/rat) *Colorimetric*	ELISA	Colorimetric
AS-55153-H	Sensolyte® Anti-Human MOG (1-125) Human IgG Specific ELISA Kit *Colorimetric*	ELISA	Colorimetric
AS-55153-M	Sensolyte® Anti-Human MOG (1-125) Mouse IgG Specific ELISA Kit *Colorimetric*	ELISA	Colorimetric

Catalog#	Kit	Assay type	Detection
AS-55156	Sensolyte® Anti-Mouse MOG (1-125) IgG Quantitative ELISA Kit *Colorimetric*	ELISA	Colorimetric
AS-55157	Sensolyte® Anti-Rat MOG (1-125) IgG Quantitative ELISA Kit *Colorimetric*	ELISA	Colorimetric
AS-55523	Sensolyte® Anti-PLP (178-191) IgG Quantitative ELISA Kit (mouse) *Colorimetric*	ELISA	Colorimetric
AS-55524	Sensolyte® Anti-PLP (139-151) IgG Quantitative ELISA Kit (mouse) *Colorimetric*	ELISA	Colorimetric

Recombinant Proteins

Catalog#	Recombinant Protein	Size
AS-55150-100	Recombinant mouse MOG (1-125)	100 µg
AS-55150-500	Recombinant mouse MOG (1-125)	500 µg
AS-55150-1000	Recombinant mouse MOG (1-125)	1000 µg
AS-55152-100	Recombinant rat MOG (1-125)	100 µg
AS-55152-500	Recombinant rat MOG (1-125)	500 µg

Catalog#	Recombinant Protein	Size
AS-55152-1000	Recombinant rat MOG (1-125)	1000 µg
AS-55158-100	Recombinant human MOG (1-125)	100 µg
AS-55158-500	Recombinant human MOG (1-125)	500 µg
AS-55158-1000	Recombinant human MOG (1-125)	1000 µg

Peptides

Catalog#	Peptide	Size
AS-22916	MBP (4-14), bovine, MBP (3-13), mouse	5 mg
AS-62081	MBP (69-85), guinea pig	1 mg
AS-29211	MBP (84-96), mouse, MBP (86-98), guinea pig, MBP (87-99), human	1 mg
AS-62731	MBP (85-98), guinea pig, MBP (86-99), human, MBP (83-96) mouse	1 mg
AS-62739	MBP (85-106) guinea pig, MBP (86-107), human	1 mg
AS-60977	MBP (88-104), guinea pig, MBP (89-105), human, MBP (86-102), mouse	1 mg

Catalog#	Peptide	Size
AS-28231	Biotin-LC-MBP Derivatized Peptide	1 mg
AS-65470	Peripheral Myelin Protein P0 (180-199), mouse	1 mg
AS-65472	Peripheral Myelin Protein P2 (53-78), bovine	1 mg
AS-63918	MOG (35-55), human	1 mg
AS-60130-1	MOG (35-55), mouse, rat	1 mg
AS-63912	PLP (139-151)	1 mg
AS-62741	PLP (178-191)	1 mg



Related Products in Multiple Pathways

Assay kits

Catalog#	Kit	Assay type	Detection
AS-72226	Sensolyte® 520 ADAM10 Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72244	Sensolyte® Transglutaminase Activity Assay Kit *Colorimetric*	Enzymatic Activity	Colorimetric
AS-72245	Sensolyte® Red Transglutaminase Assay Kit *Fluorimetric*	Enzymatic Activity	Fluorescent
AS-72254	Sensolyte® 520 Meprin β Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72149	Sensolyte® 520 Calpain Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72150	Sensolyte® AMC Calpain Activity Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72164	Sensolyte® 520 Cathepsin B Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent
AS-72165	Sensolyte® 440 Cathepsin B Assay Kit *Fluorimetric*	FRET-based Assay	Fluorescent

[Explore our full neuroscience product portfolio on our website](#)