

Proteinase-Activated Receptors

Key References

- Andrade-Gordon, P. et al. "Design, synthesis, and biological characterization of a peptide-mimetic antagonist for a tethered-ligand receptor." *Proc. Natl. Acad. Sci. USA* **96**, 12257-12262 (1999).
- Bernatowicz, M.S. et al. "Development of potent thrombin receptor antagonist peptides." *J. Med. Chem.* **25**, 4879-4887 (1996).
- Coughlin, S.R. "Thrombin signalling and protease-activated receptors." *Nature* **407**, 258-264 (2000).
- Dery, O. et al. "Proteinase-activated receptors: Novel mechanisms of signaling by serine proteinases." *Am. J. Physiol.* **274**, C1429-C1452 (1998).
- Ishihara, H. et al. "Protease-activated receptor 3 is a second thrombin receptor in humans." *Nature* **386**, 502-506 (1997).
- Kahn, M.L. et al. "A dual thrombin receptor system for platelet activation." *Nature* **394**, 690-694 (1999).
- Nakanishi-Matsui, M. et al. "PAR3 is a cofactor for PAR4 activation by thrombin." *Nature* **404**, 609-613 (2000).
- Nystedt, S. et al. "Molecular cloning of a potential proteinase activated receptor." *Proc. Natl. Acad. Sci. USA* **91**, 9208-9212 (1994).
- Rasmussen, U.B. et al. "cDNA cloning and expression of a hamster alpha-thrombin receptor coupled to Ca²⁺ mobilization." *FEBS Lett.* **288**, 123-128 (1991).
- Vergnolle, N. et al. "Proteinase-activated receptors in inflammation, neuronal signaling and pain." *Trends Pharmacol. Sci.* **22**, 146-152 (2001).
- Vu, T.K. et al. "Molecular cloning of a functional thrombin receptor reveals a novel proteolytic mechanism of receptor activation." *Cell* **64**, 1057-1068 (1991).
- Xu, W.-F. et al. "Cloning and characterization of human protease-activated receptor 4." *Proc. Natl. Acad. Sci. USA* **95**, 6642-6646 (1998).

Overview

The search for a functional thrombin receptor, using expression cloning methods, led to the discovery of a G protein-coupled receptor that mediates the cellular actions of thrombin on platelets and endothelial cells. The distinguishing property of the thrombin receptor, compared with other G protein-coupled receptors, relates to its lack of a circulating agonist and its unique mechanism of stimulation that involves the proteolytic unmasking of an activating N-terminal sequence buried in the "pro-receptor". The "revealed" N-terminal sequence acts as a "tethered" or anchored receptor-stimulating ligand. Remarkably, it was discovered that the Proteinase Activated Receptor for thrombin (now termed, PAR₁) can be activated in the absence of thrombin by relatively short peptides based on the sequence of the proteolytically revealed N-terminal domain (so-called Thrombin Receptor-Activating Polypeptides, or TRAPs). The TRAPs (now termed, PAR₁-Activating Peptides, or PAR₁APs) were used as surrogate activators of PAR₁ to evaluate the potential actions of thrombin in tissues wherein the effects of the proteinase itself might be difficult to interpret. The TRAPs also served as a basis for the development of PAR₁-targeted antagonists.

Results with the originally designed PAR₁-APs, including peptide structure-activity studies, interspecies studies of platelet aggregation and the development of PAR₁-knockout mice, clearly pointed to the existence of other members of the PAR-family. Thus, the serendipitous discovery of a second G protein-coupled proteinase-activated receptor that could be stimulated by trypsin, but not thrombin (PAR₂), was not entirely unexpected. In addition to

having considerable homology with PAR₁, PAR₂ bore the hallmark of PAR₁ in terms of the tethered-ligand mechanism for its activation by trypsin. Like PAR₁, PAR₂ can be activated by short peptides (e.g. SLIGRL... from rat PAR₂) based on the N-terminal "tethered-ligand" sequence revealed by trypsin. Whereas the PAR₂AP, SLIGRL-NH₂ cannot activate PAR₁, it was quickly realized that the originally described PAR₁APs (or TRAPs) could activate PAR₂; and that peptide structure-activity studies were required to design PAR₁APs that could selectively activate PAR₁, without activating PAR₂. Because of the cross-reactivity of PAR₁-targeted ligands with PAR₂, receptor-selective antagonists, binding probes have proved difficult to synthesize. For instance, the peptide PAR₁ antagonists that can block thrombin-mediated platelet activation have been shown to be agonists for PAR₂. Nonetheless, a non-peptide PAR₁ antagonist (RWJ 56110) has recently been described that can block both thrombin- and peptide-mediated receptor activation.

The discovery of PAR₂ did not, however, explain the activity of thrombin in murine platelets that lack PAR₁. The absence of PAR₁ in murine platelets prompted a continued search for more thrombin receptors, resulting in the discovery of two more family members, PAR₃ and PAR₄, each of which has a unique thrombin-revealed tethered ligand. Unexpectedly, PAR₃ does not appear to signal itself, but rather acts as a cofactor for the activation of PAR₄. The physiological role for PAR₃ is at present somewhat of an enigma. Synthetic peptides modeled on the thrombin-revealed sequence of PAR₃ (e.g. TFRGAP...) do not activate PAR₃, but very

likely would activate both PAR₁ and PAR₂. The PAR₄ tethered ligands (murine-GYPGKF; human-GYPGQV) fail to activate other PARs, but are of limited value as agonists for physiological studies because of their low potencies (active in the 100-400 μM range).

In summary, four proteinase-activated receptors that are regulated by a proteolytically-revealed tethered ligand mechanism are known at present. Apart from the recognized roles for PARs 1, 3 and 4 in regulating platelet and endothelial cell function, the precise physiological functions for these novel receptor systems have yet to be determined. Preliminary data point to prominent roles for the PARs in physiological processes ranging from inflammation to pain sensation and regulation of the vascular, pulmonary and gastrointestinal systems.

Proteinase-Activated Receptors

CURRENTLY ACCEPTED NAME	PAR ₁	PAR ₂	PAR ₃	PAR ₄
ALTERNATE NAMES	Thrombin receptor PAR-1	PAR-2	Thrombin receptor PAR-3	Thrombin receptor PAR-4
STRUCTURAL INFORMATION	425 aa (human)	397 aa (human)	374 aa (human)	385 aa (human)
AGONIST PROTEINASES	Thrombin (T 7513 (b), T 1063 (h)) > Trypsin (T 1426) >> Plasmin (P 4895)	Trypsin (T 1426) ^a Tryptase (T 7063) Trypsin-2 Matriptase/MT-serine protease 1 Factor Xa (F 2027)/TF/Factor VIIa	Thrombin (T 7513 (b), T 1063 (h)) >> Trypsin (T 1426) > Factor Xa (F 2027)	Thrombin (T 7513 (b), T 1063 (h)) ≈ Trypsin (T 1426) Cathepsin G (C 4428); Factors VIIa/X (F 4003)
SUBTYPE SELECTIVE AGONISTS	TFRIFD ^b TFLLR-NH ₂	SLIGRL-NH ₂	Thrombin cleaves, but does not activate	GYPGKF-NH ₂ ^c GYPGQV-NH ₂ AYPGKF-NH ₂
PUTATIVE ANTAGONISTS	BMS 200261 ^d Mercaptopropionyl-Phe-Cha-Arg-Lys- Pro-Lys-Pro-Asn-Asp-Lys-NH ₂ ^e RWJ 56110			
SIGNAL TRANSDUCTION MECHANISMS	G _{q/11} (increase IP ₃ /DAG) G _i (cAMP modulation) G _{12/13} (actin rearrangement)	G _{q/11} (increase IP ₃ /DAG) G _i (cAMP modulation)		G _{q/11} (increase IP ₃ /DAG)

Abbreviation

BMS 200261: Transcinnamoyl-p-fluoroPhe-p-guanidinoPhe-Leu-Arg-NH₂

RWJ 56110: N-((1S)-3-Amino-1-[[[(phenylmethyl)amino]carbonyl]propyl]-a-[[[1-(2,6-dichlorophenyl)methyl]-3-(1-pyrrolidinylmethyl)-1H-indol-6-yl]amino]carbonyl]amino)-3, -difluoro-benzenepropanamide

TF: Tissue factor

b: bovine

h: human

FOOTNOTES

a Thrombin is inactive.

b TFRIFD is the *Xenopus* thrombin receptor tethered-ligand domain (the human PAR₁ tethered-ligand domain sequence SFLLRN activates both PAR₁ and PAR₂ receptors).

c AYPGKF is 10-fold more potent than the natural human (GYPGQV) or murine (GYPGKF) tethered ligand sequences, which are also PAR₄ selective.

d May also act at PAR₂.

e Antagonizes PAR₁; acts as an agonist at PAR₂.