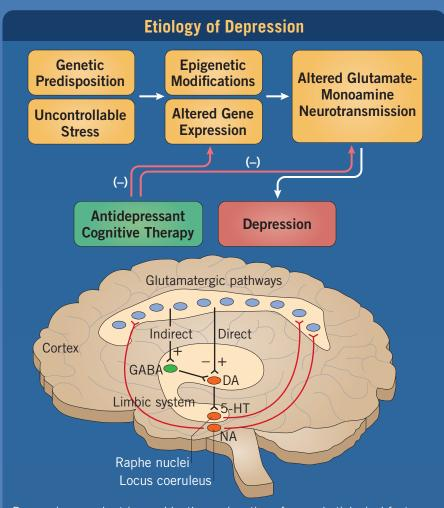
Antidepressants: Current and Future Targets

Phil Skolnick¹, Piotr Popik^{2,3} and Ramon Trullas^{4,5}

¹National Institute on Drug Abuse, National Institutes of Health, Bethesda, MD 20892 ²Institute of Pharmacology, Polish Academy of Sciences, Smetna 12, 31-343 Kraków, Poland ³Faculty of Health Sciences, Medical College, Jagiellonian University, Kraków, Poland

⁴Neurobiology Unit, Institut d'Investigacions Biomèdiques de Barcelona, Consejo Superior de Investigaciones Científicas, Institut d'Investigacions Biomèdiques August Pi i Sunyer, Rosselló 161, 08036 Barcelona, Spain ⁵Centro de Investigación Biomédica en Red sobre Enfermedades Neurodegenerativas (CIBERNED), Spain

Major depressive disorder (MDD), often referred to as major depression, is characterized by the core symptoms of depressed mood and a loss of interest and/or pleasure. Other symptoms that may be manifested include significant weight changes (loss or gain), sleep disturbances (insomnia or hypersomnia), fatigue, diminished ability to think or concentrate, feelings of worthlessness or guilt, recurrent thoughts of death or suicide, and psychomotor agitation or retardation. In addition to producing clinically significant distress, a major depressive episode is almost uniformly accompanied. by some degree of social and/or occupational impairment, negatively impacting quality of life and contributing to the societal burden associated with loss of work and health care costs. In developed countries approximately 15% of the population has been affected by MDD in their lifetime and around 5% has suffered a major depressive episode during the last year^[1]. The risk of developing MDD is almost twice as high in women^[2] and the majority of people affected with MDD do not receive standard treatment^[3].



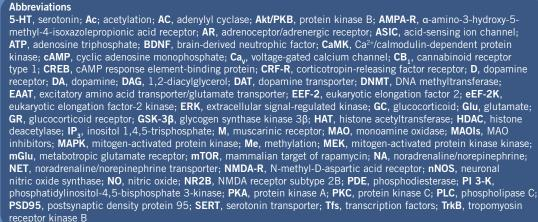
Depression may be triggered by the conjunction of several etiological factors which are capable of altering the function of brain areas associated with mood, such as the limbic system. There appears to be a strong genetic contribution to depression, and although a number of gene candidates have been reported, none have been irrefutably identified. Stressful life experiences may converge to modify gene expression that results in enduring changes in neurotransmission in limbic areas of the brain. Antidepressant drugs and cognitive therapy limit the negative effects of these etiological factors.

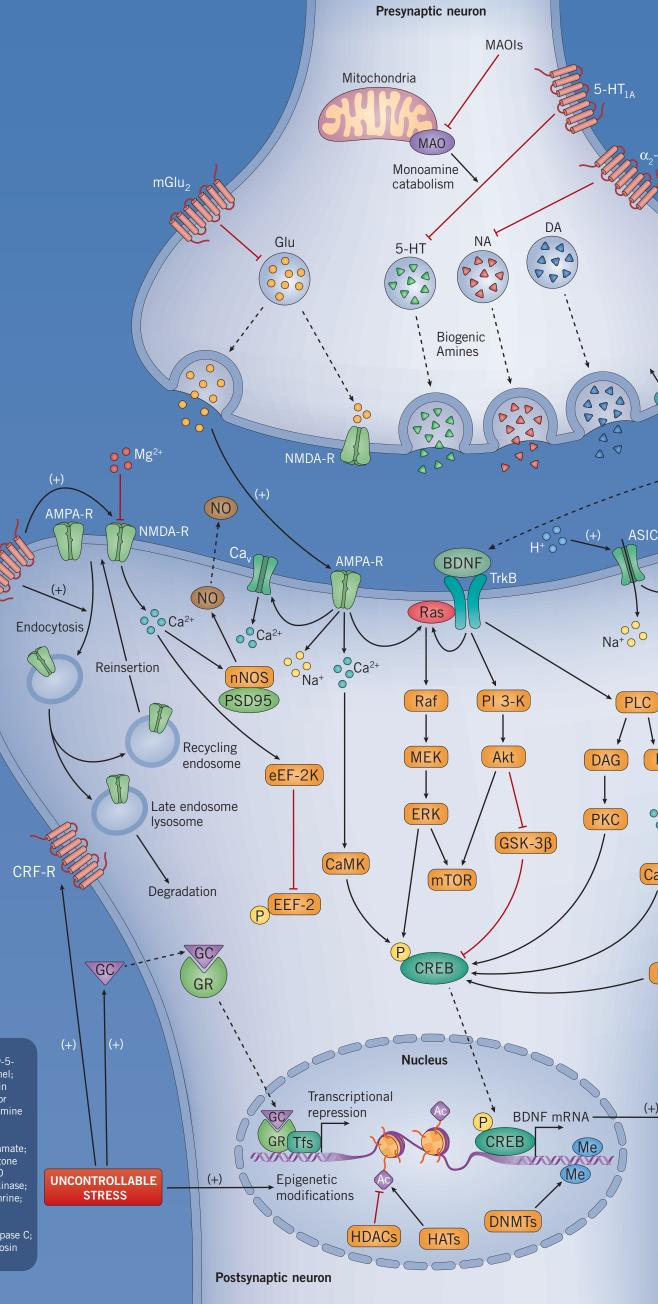
mGlu

Pharmacology of Antidepressant Drugs

All currently approved antidepressants (ADs) increase the synaptic availability of one or more of the biogenic amine transmitters: noradrenaline, dopamine, and serotonin. These ADs act either by inhibition of monoamine oxidase (MAO), an enzyme responsible for the metabolism of biogenic amines, or by inhibiting one or more of the transport proteins that remove these biogenic amines from the synaptic cleft. The latter mechanism is the principal means of terminating the actions of these neurotransmitters. The majority of ADs in current use selectively block the uptake of serotonin and/or noradrenaline. In well controlled, double blind clinical trials, biogenic amine-based ADs require \geq 2 weeks of treatment before providing meaningful relief. This "therapeutic lag" has been hypothesized to represent neuroplastic changes that must precede a reduction in depressive symptomatology.

The objective of identifying novel targets that bypass the aminergic synapse is to develop drugs that are both more effective and rapid acting than currently approved agents. Clinical studies performed over the past decade with NMDA antagonists like ketamine and traxoprodil^[4], and the muscarinic antagonist scopolamine^[5], have demonstrated that it is possible to produce a rapid and robust therapeutic response in depressed individuals, many of whom do not respond to biogenic amine based agents. While many promising targets have been identified, a high safety bar will be demanded of new therapeutics as long as alternative therapies, however imperfect, are available.







TOCRIS a biotechne brand

Neurochemicals | Signal Transduction Agents | Peptides | Biochemicals

www.tocris.com

Astrocyte EAAT M₁ Receptor ASIC1a Ca²⁺ IP₃ cAMP) ATP PDE

Targets for the Development of Antidepressants

Presynaptic targets:

- Receptors that regulate neurotransmitter release such as the 5-HT₁₄, α₂ adrenoceptor, mGlu₂ and NMDA receptor.
- MAO, an enzyme that catabolizes monoamines.
- Neurotransmitter transporters, exemplified by biogenic amine transporters. Inhibiting these transporters increases the availability of neurotransmitter at the synaptic cleft, a key action in the mechanism of most currently prescribed antidepressants.

Postsynaptic targets:

Targets for antidepressant action include both ionotropic and metabotropic receptors, as well as the signal transduction pathways activated by these receptors. These pathways converge in regulating gene expression, release of neurotrophic factors, and synaptic plasticity that may contribute to the mechanism of antidepressant action.

mGlu_{1/5} receptors regulate NMDA receptor function by modulating its trafficking and cell surface expression. The NMDA receptor is blocked by Mg²⁺ at resting membrane potential when the membrane is depolarized, this block is relieved, allowing Ca^{2+} to enter the cell. Ca^{2} participates in a number of cellular cascades, including nitric oxide synthesis, which functions as a retrograde neuronal messenger. Ca²⁺ can also enter the cell via voltage-dependent Ca²⁺ channels (Ca.) where it activates CaMK. Ca. function can be modulated by other ion channels sensitive to glutamate such as the AMPA receptor. Both AMPA receptors and the TrkB receptor (the receptor for BDNF) influence the activity of signal transduction pathways, such as the Ras-Raf-MEK-MAPK and PI 3-K-Akt-GSK-3β signaling pathways, which may be involved in the mechanism of action of certain antidepressant drugs.

Other metabotropic receptors such as 5-HT₂, a, adrenoceptors and M, muscarinic receptors, stimulate PLC and second messengers through $G_{q/11}$ signaling. G_s and G_i proteins are activated by β -adrenoceptors, $D_{1,2,5}$ receptors and 5-HT_{1A,4,5,7} receptors. These G proteins modulate AC activity, which regulates the level of cAMP and thus PKA activity. cAMP concentrations are also affected by drugs that influence PDE enzymatic activity. Scopolamine is a muscarinic receptor antagonist that inhibits M₁-M₅ with similar affinity; studies with both knock-out mice and selective pharmacological agents ^[6] indicate that the M, and to a lesser extent M_a receptor subtypes may be responsible for the rapid and robust antidepressant effects of this drug.

Many of these signal transduction pathways (CaMK, ERK, GSK-3β) appear to converge in phosphorylating transcription factors such as CREB, a transcription factor that binds to DNA sequences (cAMP response elements (CRE)), before inducing transcription and the expression of certain genes in the nucleus of the neuron. Such changes in gene expression induced by antidepressants can modify the genetic and epigenetic alterations caused by inescapable life experiences, as well as other genetic factors that are believed to be involved in the pathophysiology of depression. For example the rapid acting antidepressants (ketamine scopolamine) have been associated with rapid changes in synaptic morphology, including increases in spine density and spine head diameter

Compounds clinically validated or with preclinical evidence:

- Acetylcholine muscarinic receptor antagonists: **scopolamine**
- NMDA ion channel blockers: AZD 6765, ketamine, Mg²⁺, memantine, Zn²⁺
- NMDA receptor (glycine site) partial agonists: ACPC, **D-cycloserine**, **Glyx-13**, HA-966
- NMDA receptor (glycine site) antagonists: 4-chlorokynurenine. 5.7-dichlorokynurenic
- acid 1-701.324 • Selective NR2B antagonists: CP 101, 606 (traxoprodil), eliprodil, ifenprodil, MRK-0657
- Compounds highlighted in bold are clinically validated

Products available from Tocris acting on antidepressant pathways

PKA (+) BDNFT	Monoamine Oxidase	5-HT₄: BIMU 8, GR 113808, GR 125487,	Vasopressin Receptors
	Lazabemide, Moclobemide, Pirlindole	RS 67333	[Arg8]-Vasopressin, SR 49059
	Glutamate Receptors	5-HT₅: SB 699551	Tachykinin Receptors
	NMDA: D-AP5, CPP, 5,7-Dichlorokynurenic acid,	5-HT ₇ : AS 19, SB 258719, SB 269970	GR 73632, L-733,060, RP 67580
	DQP 1105, Ifenprodil, Ketamine, L-689,560,	Dopamine Receptors	Signal Transduction
	Memantine, (+)-MK 801, Ro 25-6981	D ₁ /D ₅ : SCH 23390, SCH 39166, SKF 81297	AC: Forskolin, NKH 477, SQ 22536
	AMPA: (S)-AMPA, Aniracetam, CX 546,	D ₂ : L-741,626, (-)-Quinpirole, Sumanirole	Akt: API-1, 10-DEBC, GSK 690693
	Cyclothiazide, GYKI 53655, Naspm, S 18986	Neurotransmitter Transporters	Ca ²⁺ Signaling: A23187, Ionomycin
	mGlu _{1/5} : FTIDC, JNJ 16259685, LSN 2463359,	SERT: Citalopram, Fluoxetine, Sertraline	CAMK: A 484954, KN 93, STO-609
	LY 367385, MPEP, MTEP, Ro 67-7476	DAT: Bupropion, GBR 12909	DNMTs: Decitabine, SGI 1027, Zebularine
	mGlu ₂ : LY 341495, LY 354740, LY 379268, LY 487379	NET: Reboxetine, Tomoxetine	GSK-3: BIO, SB 216763, SB 415286
	Glutamate Release: Lamotrigine, Riluzole	EAAT: Dihydrokainic, DL-TBOA, TFB-TBOA	HDACs: FK 228, SAHA, Trichostatin A
	Muscarinic Receptors	Ion Channels	MEK: BIX 02189, PD 0325901, PD 98059,
	Scopolamine, VU 0255035, Xanomeline	Ca _v : ω-Conotoxin GVIA, Mibefradil, SNX 482	SL 327, U0126
	Adrenergic Receptors	ASIC: Psalmotoxin 1	mTOR: KU 0063794, Rapamycin, Temsirolimus,
	a₁-AR: A 61603, Cirazoline, (<i>R</i>)-(-)-Phenylephrin, Prazosin	TASK-3: ML 365	Torin 1
		TREK: BL 1249, Spadin	NO: N ^ω -Propyl-L-arginine, SNAP
	α ₂ -AR: Dexmedetomidine, Imiloxan, RS 79948	GR	PDE: IBMX, Rolipram, Sildenafil
	β-AR: Isoproterenol, Propranolol, SR 58611A	Fluticasone, Mifepristone	PI 3-K: LY 294002, Wortmannin, 740 Y-P
	TrkB Receptor	CRF-R	PKA: H 89, KT 5720, PKI 14-22
	ANA 12, BDNF, LM 22A4	Astressin 2B, CP 376395, CRF (human, rat),	PLC: D609, U 73122
	Serotonin Receptors	NBI 35965, Stressin I	Raf: GW 5074, SB 590885
	5-HT _{1a} : 8-Hydroxy-DPAT, Xaliproden	FGFR	
	5-HT ₂ : Clozapine, MDL 100907, TCB-2	PD 173074, SU 5402, SU 6668	
	Deferences		Michalitation Destant Transfer Discourse (Coli 0000
	References:	at al Appy Roy Clip Revokal 2007	[4] Skolnick, P., et al. Trends Pharmacol.Sci. 2009
		et al. Annu.Rev.Clin.Psychol. 2007	[5] Drevets, W.C., <i>et al. Biol.Psychiatry</i> 2013

[2] Compton, W.M., et al. Am.J.Psychiatry. 2006[6] Witkin, J.M., et al. J.Pharmacol.Exper.[3] Kessler, R.C. and Bromet, E.J., Annu.Rev.Public Health. 2013[7] Voleti, B., et al. Biol.Psychiatry 2013