

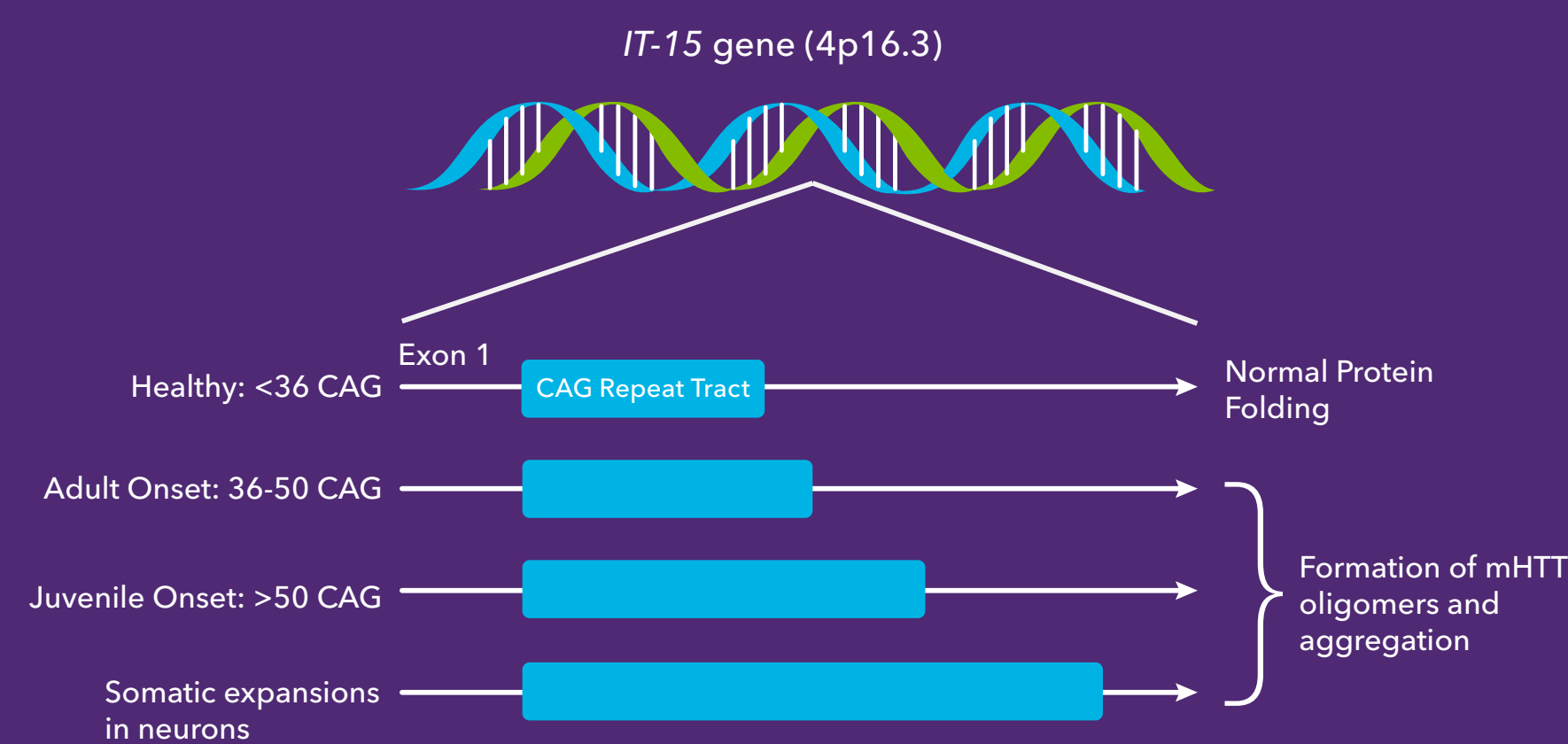
Huntington's Disease: Pathophysiology and Clinical Prospects

Nina Stöberl and Nicholas D. Allen, School of Biosciences, Cardiff University, The Sir Martin Evans Building, Museum Avenue, Cardiff, UK, CF10 3AX

Huntington's disease (HD) is a severe monogenic neurodegenerative disorder. It is caused by a dominantly inherited CAG trinucleotide repeat expansion in the huntingtin (HTT) gene, which encodes for an elongated glutamine stretch in the HTT protein. The progressive brain degeneration in HD is characterized by the prevalent loss of GABAergic medium spiny neurons (MSN) in the striatum¹. Clinical features include progressive motor dysfunctions and cognitive impairments².

Genetic and Clinical Manifestations

The monogenetic defect responsible for HD is a CAG trinucleotide repeat expansion in exon 1 of the IT15/HTT gene on chromosome 4, encoding an expanded polyglutamine (PolyQ) repeat in the huntingtin protein, which forms mutant HTT (mHTT). Individuals with less than 36 CAG repeats remain unaffected, however the presence of more than 40 CAG repeats consistently causes disease within a normal lifespan³. The number of CAG repeats inversely correlates with the age of disease onset as well as progression of motor and cognitive deficits. Clinical features typically manifest around the fourth decade. The most common motor behavior symptom of HD is chorea, which is characterized by unwanted rapid and short-lasting movements in the limbs, face, and trunk. In addition, patients exhibit severe cognitive and psychological symptoms which worsen with time².

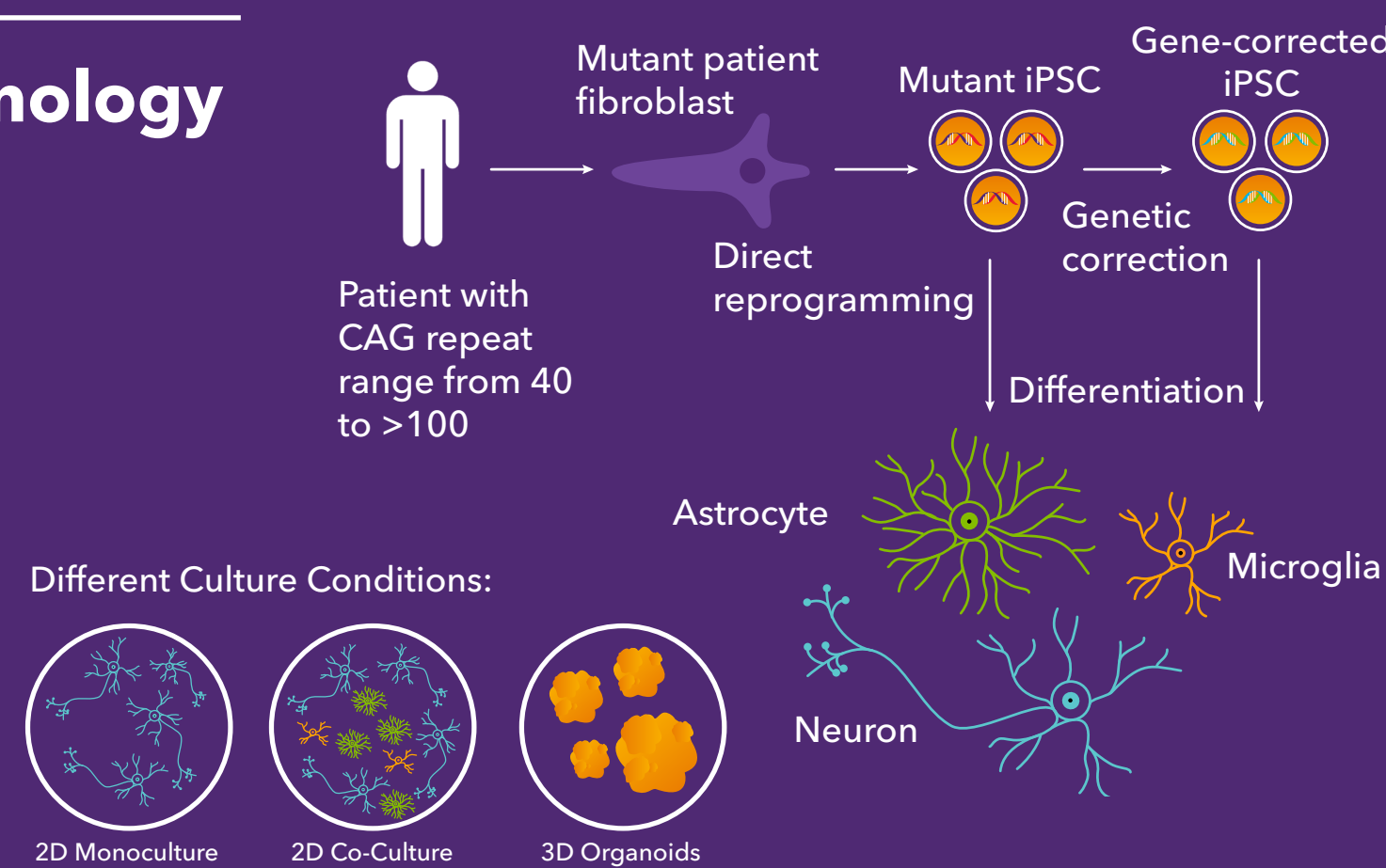


Focus on DNA Repair

Following inheritance of an expanded CAG repeat, a process of repeat expansion continues in somatic cells throughout life; higher repeat expansion rates correlate with earlier age of onset and more severe disease progression⁴. Genome-wide association studies (GWAS) have identified genes involved in DNA damage and mismatch repair pathways as genetic modifiers of disease progression and are now directly implicated in the CAG repeat expansion process⁵. This provides an exciting opportunity for HD treatment, as drugs that target identified DNA repair mechanisms have the potential to slow or even reverse somatic CAG expansion, thereby providing a therapeutic delay to disease onset^{5,6}.

HD Modeling with iPSC Technology

Somatic cells can be reprogrammed to a pluripotent state by the exogenous introduction of pluripotency associated transcription factors such as OCT4, KLF4, SOX2 and c-MYC¹³. The generated human induced pluripotent stem cells (iPSCs) can differentiate into cells of all germ layers, including brain cells like neurons, astrocytes and microglia. iPSCs can be derived from HD patients in order to retain the patient's genetic background and are accessible for gene editing, e.g. to correct the mHTT to produce isogenic control cells¹⁴. Nowadays, iPSC-derived cells can be studied in different culture conditions, ranging from 2D monocultures to 3D brain-like organoids¹⁵.



HTT Functions

The HTT protein has several functions within the cell. HTT interacts with various organelles and is involved in mediating:

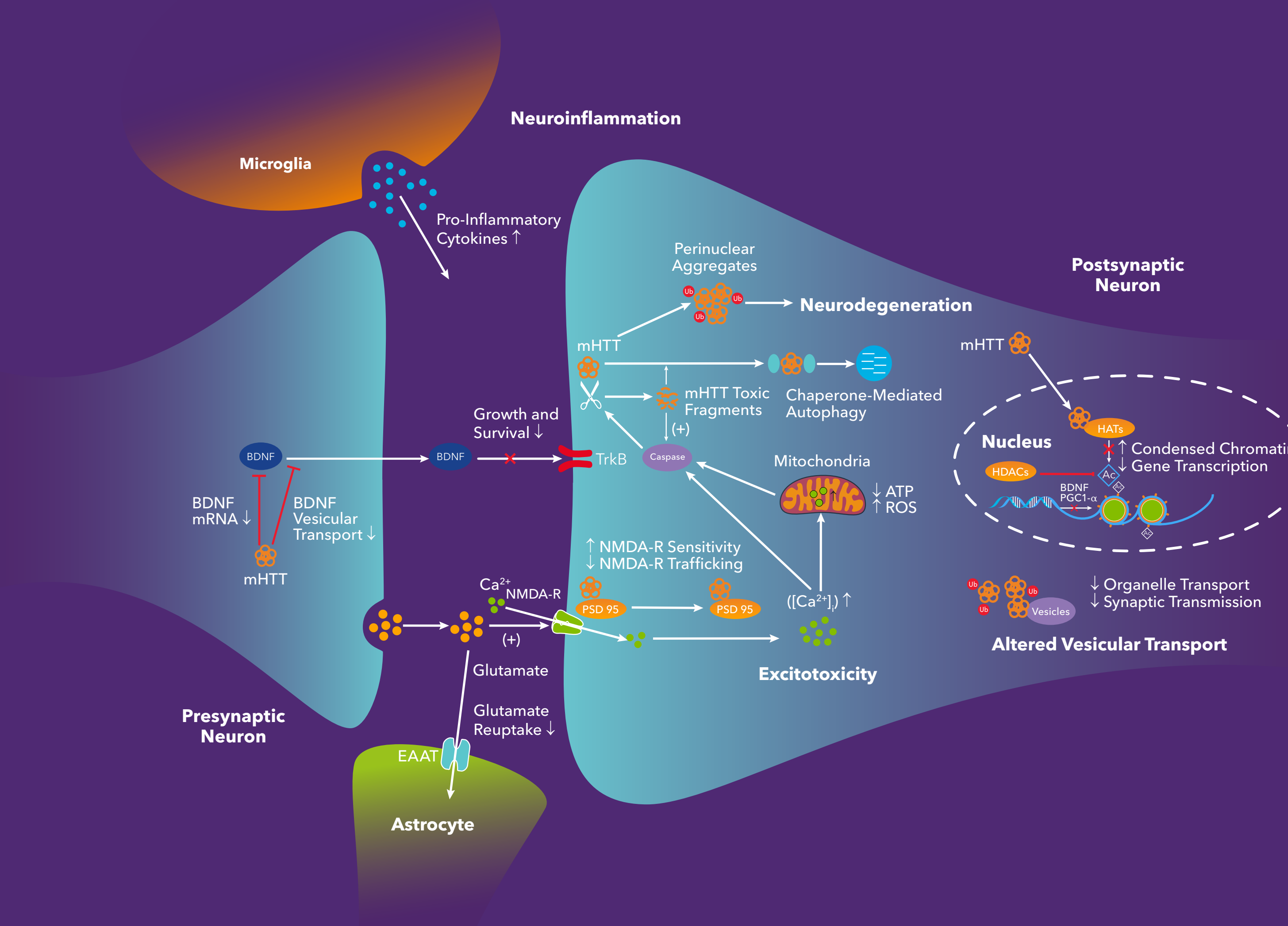
- Chromatin remodeling and gene transcription^{7,8}
- Neuroprotection (BDNF transcription and trafficking)^{9,10}
- Cell survival (HTT is anti-apoptotic)⁷
- Calcium homeostasis¹¹
- Mitochondrial function and trafficking¹⁰
- Intracellular trafficking and synaptic vesicle recycling^{7,8}

The reduction of HTT and coexistence of mHTT results in the disruption of these cellular processes, culminating in cell death⁸.

Therapeutic Strategies

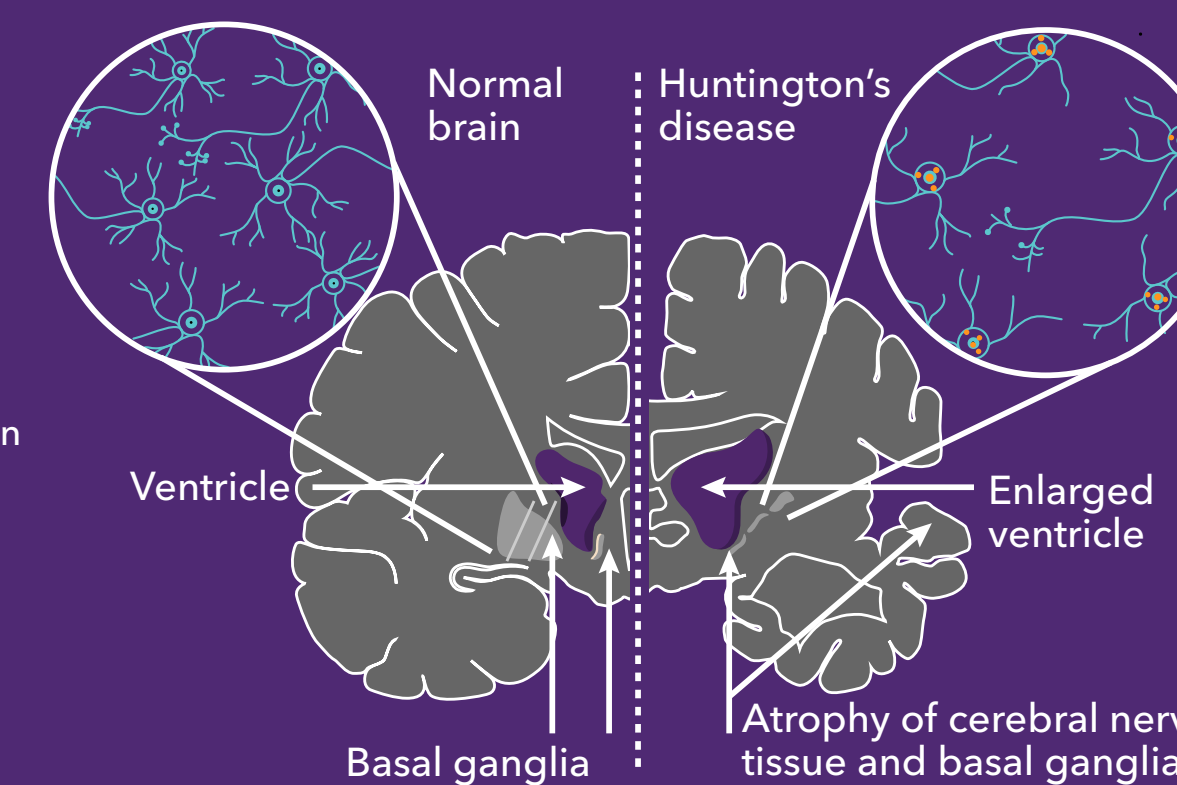
Currently, there are no treatment options for HD that can halt or reverse the disease. Commonly used symptomatic treatments attempt to improve the quality of life of affected patients by addressing the movement, cognitive and psychiatric symptoms. Novel therapeutics with the intent of slowing disease progression are being developed. These approaches target pathways specific to HD biology. Among these approaches, mHTT protein lowering therapies hold great promise¹².

Molecular Events in Neurons



Huntington's Disease Brain Pathology

The left hemisphere shows a normal healthy brain, while the right hemisphere shows HD-specific histopathological hallmarks. The most striking feature of HD is neuronal loss in the striatum along with enlarged lateral ventricles. In progressed stages of the disease, degeneration also occurs in the cerebral cortex¹⁶. At a molecular level the mHTT protein forms intracellular aggregates in neurons of affected brain regions that appear as inclusion bodies¹⁷. Expanding polyglutamine length in the mHTT protein induces increased generation of toxic mHTT oligomers and presence of cellular inclusion bodies.



References

1. Bates et al. (2015) *Nat.Rev.Dis.Primsers.* 1 15052
2. Tabrizi et al. (2020) *Nat.Rev.Neurol.* 16 529
3. Rubinsztein et al. (1996) *Am.J.Hum.Genet.* 59 16
4. Swami et al. (2009) *Hum.Mol.Genet.* 18 3039
5. Massey & Jones (2018) *Dis.Model.Mech.* 11 031930
6. Maiuri et al. (2021) *J.Huntingtons Dis.* 10 165
7. Saudou & Humbert (2016) *Neuron* 89 910
8. Gatto et al. (2020) *Clin.Park.Relat.Disord.* 3 100056
9. Cattaneo et al. (2005) *Nat.Rev.Neuosci.* 6 919
10. Jimenez-Sanchez et al. (2017) *Cold Spring Harb.Perspect.Med* 7 a024240
11. Kolobkova et al. (2017) *Acta Naturae* 9 34
12. Dash & Mestre (2020) *Neurotherapeutics* 17 1645
13. Takahashi & Yamanaka (2006) *Cell* 126 663
14. HD iPSC Consortium (2012) *Cell Stem Cell* 11 264
15. Hong & Tae Do (2019) *Front.Bioeng Biotechnol.* 7 400
16. Sieradzian & Mann (2001) *Neuropathol.Appl.Neurol.* 27 1
17. Arrasate & Finkbeiner (2014) *Exp.Neurol.* 238 1

Tocris Products

- | | |
|---|--|
| Trk Receptors
ANA 12
BDNF (human)
7,8-Dihydroxyflavone
LM 22A4 | Caspases
Cisplatin
Z-DEVD-FMK
Z-VAD-FMK |
| NMDA Receptors
D-AP5
DL-AP5
(R)-CPP
(RS)-CPP
Ibotenic acid
Memantine (+)-MK 801 | Calpains
Acetyl-Calpastatin (184-210) (human)
Calpeptin
PD 150606 |
| Kainate Receptors
CNOX
DNQX
Kainic acid
Topiramate | Cathepsin
CA 074
E 64
E 64d |
| Glutamate Transporters
Dihydrokainic acid
DL-TBOA
LDN 212320
TFB-TBOA | IP₃ Receptors
2-APB
(-)-Xestospingon C |
| Glutamate (Metabotropic) Receptors
CDPPB
CHPG Sodium salt
(RS)-3,5-DHPG
(S)-3,5-DHPG
L-Quisqualic acid
LY 379268
(S)-MCPG
MPEP
MTEP
VU 0360172 | Mitochondrial Permeability Transition Pore
Cyclosporin A
TRO 19622 |
| Heat Shock Proteins
17-AAG
Geldanamycin
VER 155008 | Proteasome
Lactacystin
MG 132 |
| Stem Cell Differentiation
DAPT
DMH-1
Dorsomorphin
Fluoxetine
Forskolin
IBMX
Purmorphamine
SAG
SB 431542
SU 5402 | Autophagy
Bafilomycin A1
(±)-Bay K 8644
Dexamethasone
LY 294002
Rapamycin
Taxol
Thapsigargin
Torin 1
Tunicamycin |
| iPSCs
CHIR 99021
Valproic acid sodium salt | Dopamine Receptors
Dopamine
Olanzapine
Risperidone
SCH 23390
SKF 38393
Tetrabenazine |
| Histone Deacetylases
SAHA
Trichostatin A | Adenosine A_{2A} Receptors
CGS 21680
Istradefylline
SCH 58261
ZM 241385 |
| SIRT
EX 527
Nicotinamide
Resveratrol | Neuroprotective Compounds
Minocycline
N-Acetylcysteine
amide
Riluzole |
| Histone Acetyltransferases
Anacardic acid
C 646 | Caged Compounds
MNI-caged-L-glutamate
RuBi-Glutamate
RuBi-GABA |