Cell Cycle and DNA Damage Repair

Adapted from Edition 3 of the Tocris Cancer Product Guide

is synthesized by Pol δ . PCNA is a cofactor for

both DNA polymerase δ and ϵ , where it acts as

a DNA clamp, which is important in both DNA

synthesis and repair. At the end of the termination

phase, DNA ligases form a phosphodiester bond,

which joins the DNA strands together, forming

new doubled stranded DNA.

To request a copy of the Tocris Cancer Product Guide, or to view the PDF, please visit tocris.com

TOCRIS

Neurochemicals | Signal Transduction Agents | Peptides | Biochemicals www.tocris.com

Products available from Tocris

In normal cells, each stage of the cell cycle is tightly regulated. In cancer cells, many genes and proteins that influence the progression of the cell cycle are mutated or overexpressed – they become oncogenes. The proteins/molecules involved in the regulation of the cell cycle, in particular DNA replication and DNA damage are important cancer therapeutic targets.

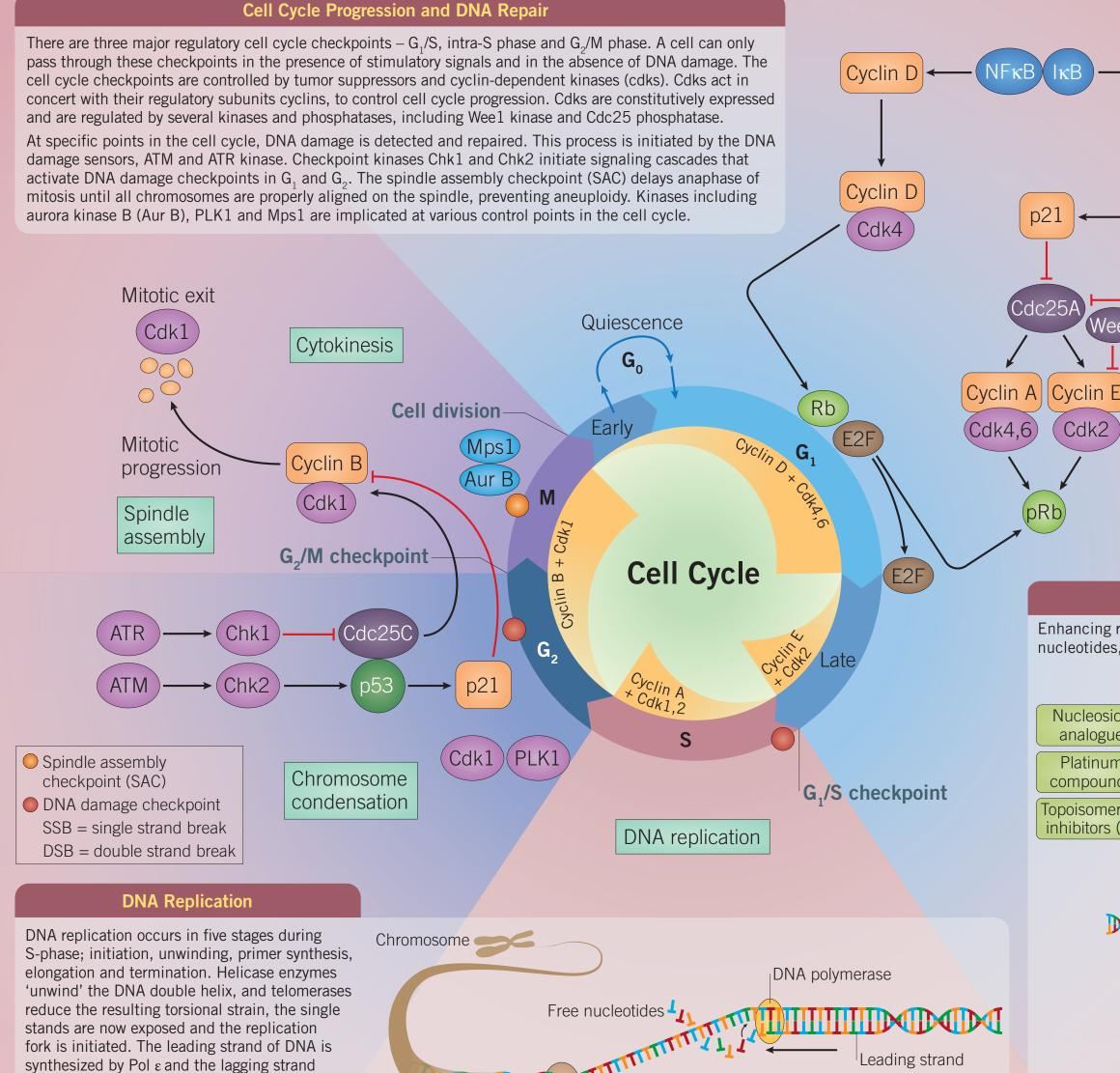
Lagging

strand

шишшиш

strand

Original (template)



Original DNA

► Thymine ► Guanine

ATM & ATR Kinase AZ 20, CGK 733, KU 55933, KU 60019, Mirin Aurora Kinases Hesperadin, TC-A 2317, VX 680, ZM 447439 Calpains Acetyl-Calpastatin (184-210) (human), Calpeptin, E 64, MDL 28170, MG 101, MG 132, PD 150606 Casein Kinase 1 D 4476, (R)-DRF053, LH 846, PF 4800567, PF 670462 Casein Kinase 2 TBB, TBCA, TMCB, TTP 22 Cdc25 Phosphatase NSC 663284, NSC 95397 Cell Cycle Inhibitors CFM 4, 10058-F4, Methotrexate, Narciclasine Checkpoint Kinases AZD 7762, CCT 241533, NSC 109555, PD 407824, PF 477736, SB 218078, TCS 2312 Cyclin-dependent Kinase Kenpaullone, PD 0332991, Purvalanol A, Purvalanol B, Ro 3306, Senexin A DNA-dependent Protein Kinase Compound 401, DMNB, NU 7026, NU 7441 DNA, RNA and Protein Synthesis 4E1RCat, L189, Mithramycin A, NSC 617145, T2AA Hsp**70** VER 155008 **Hsp90** 17-AAG IRE1 APY 29 Kinesin BRD 9876, Dimethylenastron, K 858, Monastrol, SB 743921 S-TritvI-L-cysteine Microtubules Docetaxel, Dolastatin 10, Flutax 1, Taxol, Vinblastine, Monopolar Spindle 1 Kinase AZ 3146, Mps BAY 2a, Mps1-IN-1, TC Mps1 12 p53 NSC 319726, Nutlin-3, PRIMA-1MET, RETRA, RITA Pim Kinase PIM-1 Inhibitor 2, R8-T198wt, TCS PIM-1 1, TCS PIM-1 4a Polo-like Kinase Cyclapolin 9, GW 843682X, SBE 13, TAK 960, TC-S 7005 Poly (ADP-ribose) Polymerase JW 55, NU 1025, PJ 34, XAV 939 Telomerase BIBR 1532, Costunolide, 5-TAMRA SE, TMPyP4 tosylate

ta and O'Shaughnessy (2010) Clin. Cancer Res. Lapenna and Giordano (2009) Nat. Rev. Drug Disc

Targeting Cancer Cells

Enhancing replicative stress by targeting critical DNA replication checkpoints and replication machinery, as well as depleting nucleotides, encourages fork stalling and fork collapse, which leads to mitotic catastrophe and death in cancer cells.

Proteasome

mdm2

DSBs

ATM

SSBs

DNA damage

(Chk2

Chk1

