

# Pre-profiled anti-cancer drugs

Compound	Synonym	Target
A-420983		Lck
abemaciclib	LY2835219	CDK4 / CDK6
ABT-737		BCL2 / BCL-xL / BCL-w
acalabrutinib	ACP-196	BTK
ACBI2		SMARCA2
actinomycin D	RASP-101 / Deactinomycin	transcription
adagrasib	MRTX849	KRAS G12C
adavosertib	AZD-1775	WEE1
afatinib	BIBW-2992	EGFR
alectinib	CH5424802	ALK
all-trans retinoic acid	ATRA	RAR
alpelisib	BYL719	PI3Ka
AMG-900		Aurora kinases
ART558		polymerase theta
asciminib	ABL001	ABL
AT7519		CDK
AT9283		Aurora kinases / JAK2 / JAK3
avapritinib	BLU-285	KIT / PDGFRA
avutometinib	VS-6766	MEK
axitinib	AG-013736	multikinase
AZD8055		mTOR
AZD9574		PARP1
azenosertib	ZN-c3	Wee1
batoprotafib	TNO155	SHP2
BAY 11-7085		NFKB
BAY 1217389		TTK
BAY-293		SOS1
BAY-320		BUB1
BDTX-1535	EGFR-IN-76	EGFR
belvarafenib	HM95573	pan-RAF
berzosertib	VE-822	ATR
BI 2536		PLK1
BIIB021		HSP90
binimetinib	ARRY-162	MEK
BLU-945		EGFR
BLU9931		FGFR4
bortezomib	PS-341	proteasome
BOS172722		TTK
bosutinib	SKI-606	SRC / ABL
brigatinib	AP26113	ALK / ROS1 / EGFR
brivanib	BMS-540215	multikinase
buparlisib	BKM-120	PI3K / tubulin

# Pre-profiled anti-cancer drugs

Compound	Synonym	Target
busulfan		DNA alkylating
cabozantinib	XL-184	c-MET / VEGFR
camptothecin	CPT	topoisomerase I
capivasertib	AZD5363	Pan-Akt
capmatinib	INC280	c-MET
carboplatin		DNA damage
carfilzomib	PR-171	proteasome
ceralasertib	AZD6738	ATR
ceritinib	LDK-378	ALK
cetuximab	C225	EGFR
CFI-402257		TTK
CHIR-124		CHK1
cisplatin		DNA damage
cobimetinib	GDC-0973	MEK
copanlisib	BAY 80-6946	PI3K
crenolanib	CP-868596	FLT3 / PDGFRa / PDGFRb
crizotinib	PF-02341066	ALK / ROS1 / c-MET
CYC116		Aurora kinase A / Aurora kinase B
cyclophosphamide	4-hydroxy cyclophosphamide	DNA alkylating
cytarabine		nucleoside analog
D08	2-D08	SUMOylation
dabrafenib	GSK-2118436	BRAF-monomer
dacarbazine		DNA alkylating
dacomitinib	PF299804, PF299	EGFR
dactolisib	BEZ-235	PI3K / mTOR
danusertib	PHA-739358	Aurora kinases
dasatinib	BMS-354825	SRC / ABL
daunorubicin		topoisomerase II
Debio 0123		WEE1
decitabine	NSC 127716	DNA methyltransferase
defactinib		FAK
degrasyn	WP1130	deubiquitinase
derazantinib	ARQ087/ BAL087	FGFR1 / FGFR2 / FGFR3
deruxtecan	DXd	topoisomerase I
dinaciclib	SCH-727965	CDK
docetaxel	RP56976	beta-tubulin
doxifluridine	5-dFdU	nucleoside analog
doxorubicin	Adriamycin	topoisomerase II
DT2216		BCL-xL
duvelisib	IPI-145	PI3Kd / PI3Kg
elimusertib	BAY-1895344	ATR
enasidenib	AG-221	IDH2
encorafenib	LGX818	BRAF-monomer

# Pre-profiled anti-cancer drugs

Compound	Synonym	Target
entinostat	MS-275	HDAC
entrectinib		NTRK / ROS1
epacadostat		IDO1
epirubicin	pharmorubicin	topoisomerase II
eprenetapopt	APR-246/ PRIMA-1MET	TP53
erdafitinib	JNJ-42756493	FGFR
eribulin		tubulin
erlotinib	OSI-774	EGFR
etoposide	VP-16	topoisomerase II
everolimus	RAD001	mTOR / FKBP12
exatecan		topoisomerase I
fenebrutinib	GDC-0853	BTK
fluorouracil	5-FU	nucleoside analog
fostamatinib		SYK
fulvestrant	ZD 9238	ESR
futibatinib	TAS-120	FGFR
GDC-0152		XIAP / IAP
gefitinib	ZD-1839	EGFR
gemcitabine	LY-188011	nucleoside analog
gilteritinib	ASP2215	FLT3 / ALK
GSK1070916		Aurora kinase B/ Aurora kinase C
GSK126		EZH2
GSK343		EZH2
GSK461364		PLK1
I-138		USP1-UAF1
ibrutinib	PCI-32765	BTK / EGFR
ICG-001		Wnt-pathway
idelalisib	CAL-101	PI3Kd
imatinib	STI-571	ABL
infigratinib	NVP-BGJ398	FGFR
ipatasertib	GDC-0068	pan-AKT
irinotecan	CPT-11	topoisomerase I
ivosidenib	AG-120	IDH1
JA2131		PARG
JBj-09-063		EGFR
JQ1		BET
K-975		TEAD
KSQ-4279	USP1-IN-1	USP1
KU-60019		ATM
lapatinib	GSK-572016	HER2
LB42708		farnesyltransferase

# Pre-profiled anti-cancer drugs

Compound	Synonym	Target
lenvatinib	E7080	multikinase
lifirafenib	BGB-283	RAF / EGFR
linsitinib	OSI-906	IGF1R
lirafugratinib	RLY-4008	FGFR2
lorlatinib	PF-6463922	ALK / ROS1
luminespib	NVP-AUY922	IGF1R
LY3009120	DP-4978	pan-RAF
LY3295668	AK-01	Aurora kinase A
masitinib	AB-1010	KIT / PDGFRa / PDGFRb
melphalan		DNA alkylating
mercaptopurine	6-MP	nucleoside analog
methotrexate	MTX	folate synthesis
MG-132		proteasome
midostaurin	PKC412	FLT3
miransertib	ARQ-092	pan-AKT
mirdametinib	PD0325901	MEK
mitomycin C		DNA alkylating
mitoxantrone		topoisomerase II / PKC
MK-5108	VX-689	Aurora kinase A
MK-8353	SCH900353	ERK1 / ERK2
ML-792		SAE
MLN7243		ubiquitin E1 enzyme
MLN8054		Aurora kinase A
MMAE		tubulin
mobocertinib	TAK-788, AP32788	EGFR
molibresib	I-BET-762, GSK-525762	BET
momelotinib	CYT-387	JAK1 / JAK2
MPI-0479605		TTK
Mps1-IN-1		TTK
MRK-560		gamma-secretase
MRTX1133		KRAS G12D
MS934		MEK
MYCi975	NUCC-0200975	MYC
MYCMI-6	NSC354961	MYC
naporafenib	LXH254	BRAF / CRAF
navitoclax	ABT-263	BCL2 / BCL-xL / BCL-w
nedisertib	M3814	MEK
nemtabrutinib	ARQ 531/MK-1026	BTK / TEC / SRC / MEK
neratinib	HKI-272	HER2 / EGFR
nilotinib	AMN-107	ABL

# Pre-profiled anti-cancer drugs

Compound	Synonym	Target
nintedanib	BIBF-1120	multikinase
niraparib	MK-4827	PARP1 / PARP2
NMS-P153		TTK
NMS-P715		TTK
novobiocin	NSC 2382	polymerase theta
nutlin-3a		MDM2
obatociclib	GX15-070	pan-BCL
olaparib	AZD2281	PARP1 / PARP2
osimertinib	AZD9291	EGFR
paclitaxel	NSC 125973	beta-tubulin
palbociclib	PD-0332991	CDK4 / CDK6
panobinostat	LBH589	HDAC
pazopanib	GW-786034	multikinase
PDD00017273		PARG
pemetrexed		antifolate
pemigatinib	INCB054828	FGFR
pemrametostat	GSK336595	PRMT5
pevonedistat	MLN4924	NEDD8
pexidartinib	PLX3397	CSF-1R
PHA-793887		CDK
pirtobrutinib	LOXO-305	BTK
plixorafenib	PLX8394	BRAF-monomer
ponatinib	AP-24534	ABL
pralsetinib	BLU-667	RET
prednisolone		NR3C1
quizartinib	AC-220	FLT3
rebastinib	DCC-2036	TIE2 / ABL
regorafenib	BAY-73-4506	multikinase
reversine		TTK
revumenib	SNDX-5613	menin-MLL
ribociclib	LEE011	CDK4 / CDK6
ripretinib	DCC-2618	KIT / PDGFRa
RMC-7977		RAS(ON)
rucaparib	PF-01367338	PARP1 / PARP2
ruxolitinib	INCB-18424	JAK
sacituzumab govitecan	Trodelvy, SG	TROP2
sapanisertib	MLN0128	mTOR
saruparib	AZD5305	PARP1
savolitinib		c-MET
SCH772984		ERK1 / ERK2

# Pre-profiled anti-cancer drugs

Compound	Synonym	Target
SCH900776	MK-8776	CHK1
selpercatinib	LOXO-292	RET
selumetinib	AZD-6244	MEK
SN-38	irinotecan metab.	topoisomerase I
sorafenib	BAY 43-9006	multikinase
sotorasib	AMG-510	KRAS G12C
S-trityl-L-cysteine		kinesin Eg5
sunitinib	SU-11248	multikinase
talazoparib	BMN 673	PARP1 / PARP2
tamatinib	Fostamatinib metab.	SYK
tamoxifen	ICI 46474	ER
taselisib	GDC-0032	PI3Ka / PI3Kg / PI3Kd
tazemetostat	EPZ-6438	EZH2
temsirolimus	CCI-779	mTOR / FKBP12
tepotinib	EMD-1214063	c-MET
TH588		Tubulin / MTH1
thioguanine		nucleoside analog
tivozanib	AV-951	multikinase
tomivosertib	eFT-508	MNK1 / MNK2
topotecan		topoisomerase I
tovorafenib		pan-RAF
tozasertib	VX-680 / MK-0457	Aurora kinases
trametinib	GSK-1120212	MEK
trastuzumab	Herceptin	HER2
trastuzumab deruxtecan	Enhertu, T-DXd	HER2
trastuzumab emtansine	Kadcyla, T-DM1	HER2
tucatinib	ARRY-380	HER2
ulixertinib	BVD-523	ERK1/ ERK2
UMI-77		MCL-1
UNC1999		EZH1 / EZH2
vandetanib	ZD-6474	multikinase
vecabrutinib	SNS-062	BTK
vemurafenib	PLX-4032	BRAF-monomer
venetoclax	ABT-199	BCL2
vinblastine		tubulin
vincristine	leurocristine	tubulin
vinflunine		tubulin
vociprotafib	RMC-4630	SHP2
volasertib	BI-6727	PLK1
vorinostat	SAHA / MK-0683	HDAC
XAV939		TNKS (tankyrase)
zanubrutinib	BGB-3111	BTK