

## Therapeutic approaches to enhance natural killer cell cytotoxicity: the force awakens

Richard W. Childs and Mattias Carlsten

Scientific insights into the human immune system have led to unprecedented breakthroughs in immunotherapy, and drugs and cell-based therapies that have been developed to bolster humoral and T cell immune responses represent an established and growing component of cancer therapeutics. Although NK cells have long been known to have advantages over T cells in terms of their capacity to induce antigen-independent immune responses against cancer

cells, their therapeutic potential in the clinic has been largely unexplored. Here, we present different pharmacological and genetic strategies to bolster NK cell antitumour immunity. These approaches, as well as advances in our ability to expand NK cells ex vivo and manipulate their capacity to home to the tumour, have now armed investigators with a variety of new strategies to harness the full potential of NK cell-based cancer immunotherapy in the clinic.

# Lonza

#### NK cell tumour killing

NK cells can mediate cytotoxicity through several distinct mechanisms. Degranulation is the most studied pathway, in which NK cells release cytotoxic granulae upon interaction with target cells; this is controlled by NK cell receptors such as NKG2D, DNAM1 and others, which are counterbalanced by signalling through inhibitory receptors. The 'missingself' hypothesis, formulated in the 1980s, postulated the lack of self-MHC class I molecules on target cells as a common factor leading to NK cell cytotoxicity. Subsequent research has revealed that the functional capacity of NK cells is 'tuned' by self-MHC class I molecules (known as NK cell education). In humans, KIRs and NKG2A are currently the only known receptors to mediate functional tuning. The Fc receptor CD16 can trigger potent degranulation upon interaction with antibody-coated cells (known as ADCC) without a need for simultaneous co-activation signals. Other routes by which NK cells can kill are through the death receptor pathways TRAIL-TRAILR and FAS-FASL. These receptors induce apoptosis through caspase activation inside target cells and, therefore, mediate cytotoxicity independent of both NK cell education and signalling from receptors controlling NK cell degranulation.

#### Cellular therapies and ex vivo manipulation of NK cells

Adoptive transfer of short-term, ex vivo, IL-2-activated allogeneic NK cells has been shown to induce clinical responses in patients with AML and multiple myeloma. Administration of IL-2 after adoptive cell transfer can further promote in vivo expansion of infused NK cells. However, interest has shifted to the use of IL-15 to avoid expanding T\_\_ cells. Numerous methods have also been developed to expand NK cells ex vivo, which enables the use of multiple large-number infusions of highly activated NK cells.

Genetic manipulation of NK cells before adoptive transfer may allow for the optimization of in vivo persistence, homing to tumours and tumour cytotoxicity; NK cells engineered to express an anti-CD19 CAR are in clinical trials. Genetic manipulation of primary NK cells using viral vectors is currently inefficient, but new GMP-compliant methods to reprogramme NK cells using mRNA electroporation offer rapid and cost-efficient ways to explore a wide range of genetic approaches to enhance NK cell immunotherapy. A complementary approach is the viral transduction of NK cell lines using adenoviral and lentiviral vectors, enabling stable transgene expression. However, infusions of allogeneic NK cell lines require conditioning of the patient to avoid rapid rejection of the infused cells by the host immune system.

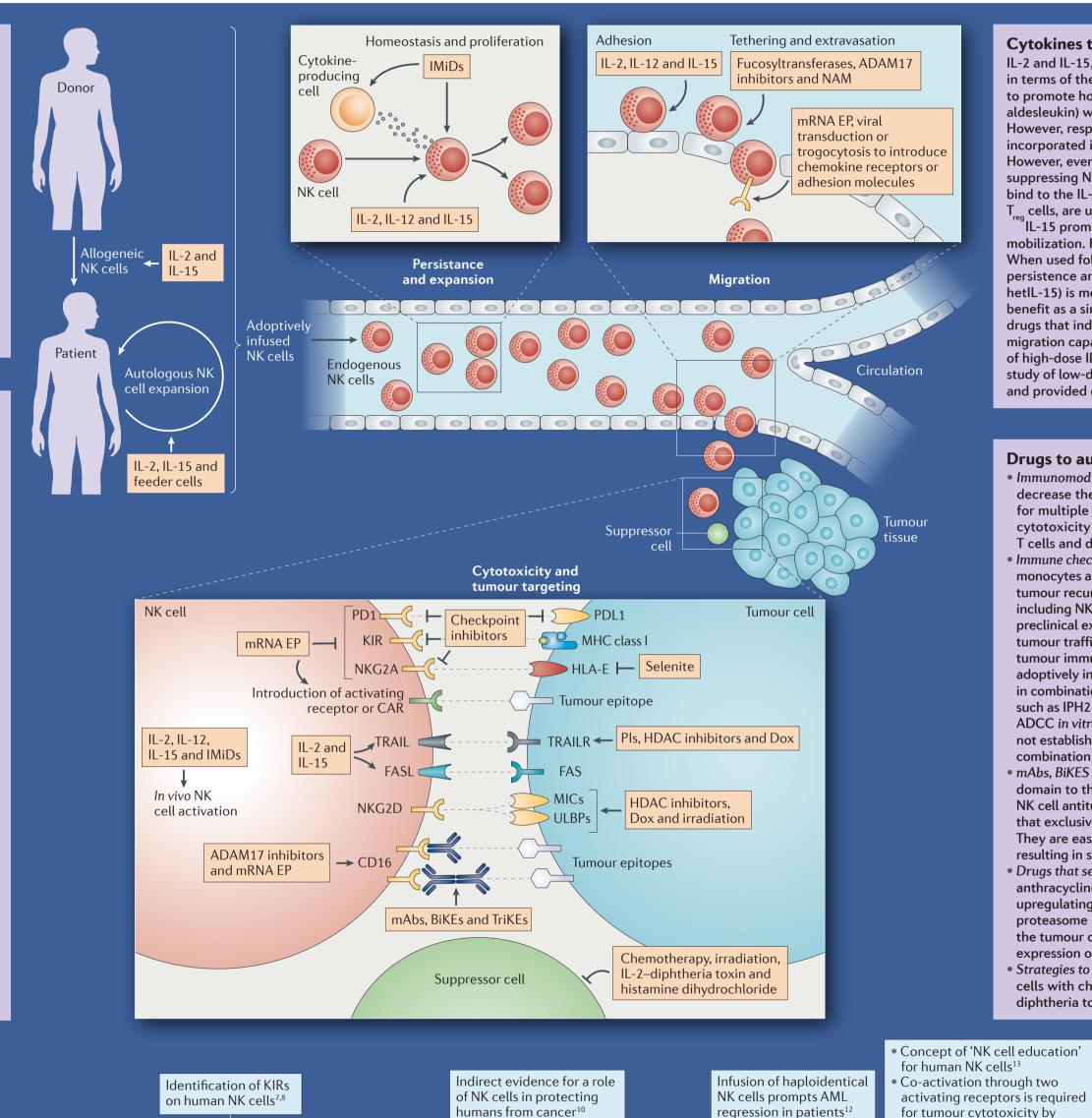
There are also several investigational strategies to manipulate ex vivo expanded NK cells with small-molecule drugs. For example, treatment with inhibitors of ADAM17 was shown to augment NK cell ADCC by preventing shedding of the CD16 receptor, and treatment of NK cells with nicotinamide enhances their expression of L-selectin, which is known to be essential for cellular trafficking.

One of the hurdles of using cytokine-activated and ex vivo expanded NK cells in patients with haematological malignancies is that NK cells express mostly non-fucosylated ligands for E-selectin, which limits their ability to home to the bone marrow. Preclinical investigations show that ex vivo fucosylation of adoptively transferred NK cells improves their antitumour effects in haematological cancers.

A novel, alternative strategy to improve NK cell migration is to alter their phenotype using mRNA EP or by culturing them on feeder cells that express homing receptors (such as CCR7), which are transferred to the NK cell membrane by trogocytosis.

NK cells shown to

mediate ADCC⁴



Cytokines to boost NK cell persistence, expansion, cytotoxicity and migration IL-2 and IL-15, as well as the pro-inflammatory cytokine IL-12, are currently being characterized in terms of their ability to stimulate NK cell antitumour immunity in humans. IL-2 has been shown to promote homeostasis, proliferation and cytotoxicity of NK cells, and rIL-2 (also known as aldesleukin) was the first cytokine used clinically to boost immune responses in cancer patients. However, responses were limited and toxicity substantial. Low-dose IL-2 therapy has been incorporated into clinical trials to support the in vivo persistence of adoptively infused NK cells. However, even ultra-low doses of IL-2 have been shown to stimulate the expansion of host T<sub>max</sub> cells, suppressing NK cell proliferation and cytotoxicity. New variants of IL-2, constructed to selectively bind to the IL-2R $\beta$  subunit expressed on NK cells rather than the IL-2R $\alpha$  subunit expressed on

IL-15 promotes NK cell development, expansion and homeostasis while avoiding T cell mobilization. However, scIL-15, the first version of IL-15 to be used in the clinic, had high toxicity. When used following adoptive NK cell infusion in patients with AML, scIL-15 supported both the persistence and proliferation of NK cells. A heterodimeric IL-15–sIL-15R $\alpha$  complex (known as hetIL-15) is more potent at stimulating NK cell proliferation. Although unlikely to be of great benefit as a single agent, it may be useful in combination with adoptive NK cell infusions or with drugs that induce NK cell tumour cytotoxicity. IL-12 upregulates adhesion molecules and the migration capacity of NK cells. Despite limited efficacy and substantial toxicity, early clinical studies of high-dose IL-12 therapy in patients with cancer showed some antitumour responses. A phase I study of low-dose subcutaneous rIL-12 in healthy individuals reported an improved toxicity profile and provided evidence that this cytokine modulates NK cell migration.

#### Drugs to augment NK cell cytotoxicity and tumour targeting

• Immunomodulatory drugs. The thalidomide derivatives lenalidomide and pomalidomide can decrease the threshold for NK cell activation. Lenalidomide represents a standard therapy for multiple myeloma and myelodysplastic syndromes, and it indirectly augments NK cell cytotoxicity and proliferation by stimulating the release of IL-2 and IFN-y from surrounding T cells and dendritic cells.

Immune checkpoint inhibitors. The checkpoint protein PD1 (expressed on activated T cells, B cells, monocytes and NK cells) and its ligand PDL1 (expressed by tumour cells) have a central role in tumour recurrence and progression, as signalling through this pathway suppresses lymphocytes, including NK cells. In vitro, blockade of PD1 on NK cells augments lysis of autologous tumour cells; preclinical experiments have shown that PD1 blockade boosts NK cell-mediated ADCC and NK-cell tumour trafficking, as well as suppressing T<sub>ma</sub> cell function. To what extent these mAbs bolster tumour immunity through NK cells, and their potential to enhance the antitumour effects of adoptively infused NK cells, remains to be investigated. PD1 inhibitors are currently in clinical trials in combination with lirilumab (also known as IPH2102), a mAb targeted against KIR. Anti-KIR mAbs such as IPH2101 and IPH2102 were shown to augment NK cell-mediated lysis of tumour cells and ADCC in vitro. However, a phase II clinical trial of IPH2101 in patients with multiple myeloma did not establish efficacy. Whether anti-KIR mAbs will show efficacy in other disease settings or in combination with other therapies remains to be determined.

• mAbs, BiKES and TriKEs: Tumour-targeting mAbs can induce ADCC through binding of their constant domain to the CD16 receptor on NK cells. However, the degree to which these mAbs mediate NK cell antitumour responses is poorly characterized. BiKEs and TriKEs are engineered molecules that exclusively act through ADCC by crosslinking epitopes on tumour cells with CD16 on NK cells. They are easier to produce than mAbs and bind to a different region of the CD16 molecule, resulting in stronger NK cell ADCC. Several BiKEs and TriKEs are in preclinical investigations.

• Drugs that sensitize tumours to NK cells. Proteasome inhibitors, such as bortezomib and the anthracycline doxorubicin, can enhance the susceptibility of tumour cells to NK cell killing by upregulating TRAILR on tumour cells and increasing the activity of caspase 8. Moreover, some proteasome inhibitors and histone deacetylase inhibitors can upregulate NKG2D ligands on the tumour cell surface, and selenite was shown to enhance NK cell killing by reducing the expression of the MHC class I molecule HLA-E on tumour cells.

Memory-like NK cells,

Description of a highly

expand large numbers

of clinical grade NK cells

efficient method to

ex vivo for adoptive

infusion in humans¹⁶

2009

previously described in mice,

now described in humans<sup>17</sup>

• Strategies to inhibit suppressor cells. NK cell activity can also be enhanced by inhibiting suppressor cells with chemotherapy, irradiation or histamine dihydrochloride, or by killing them with diphtheria toxin conjugated to IL-2.

## 1975 1977 1985 **Lonza – Your Partner in Immunotherapy Research**

Lonza is committed to supporting your immunotherapy research by providing primary cells, media and both large- and small-volume transfection solutions to help you transition from research into therapy. If you're ready for the clinic, our cell therapy services might be the answer you need for development, manufacturing and testing of your cell-based therapeutics.

#### Cells and Media

NK cells described by

Kiessling et al.<sup>2</sup> and

Hebermann et al.<sup>3</sup>

Whether you need purified natural killer cells, a mixed population of PBMCs, or unprocessed bone marrow, Lonza has the optimal cell and media solution you're looking for. If you don't see the cells that you need in our catalog, we can likely isolate them for you through our Cells on Demand Service.

#### **Transfection**

The 'missing-self'

recognition

hypothesis<sup>5</sup>

Lonza's new 4D-Nucleofector™ LV Unit can transfect up to 1x10° cells without the need for re-optimization from a smaller-scale format. With this new addition to the established 4D-Nucleofector™ System, small- and large-scale transfection applications are seamless.

1995

### Clinical Cell Therapy Services

Structures of

reported<sup>6</sup>

inhibitory NK cell

surface receptor

1989

We partner with clients to develop client-specific custom protocols to meet each unique business need, from raw material requirements and process flow to storage and distribution.

Please visit <u>www.lonza.com/immunotherapy</u> for additional information or contact scientific.support@lonza.com

#### **Abbreviations**

2000

HLA identified as predominant

ligand for the NK cell inhibitory

1998

receptor NKG2A9

ADAM17, disintegrin and metalloproteinase domain-containing protein 17; ADCC, antibody-dependent cellular cytotoxicity; ALL, acute lymphoblastic leukaemia; AML, acute myeloid leukaemia; BCL, B cell lymphoma; BiKE, bispecific killer engager; CAR, chimeric antigen receptor; CCR7, CC-chemokine receptor 7; CLL, chronic lymphatic leukaemia; CXCR3, CXC chemokine receptor type 3; DNAM1, DNAX acessory molecule 1; Dox, doxorubicin; EBV-LCL, Epstein Barr virus-lymphoblastoid cell line; EGFR, epidermal growth factor receptor; EP, electroporation; FASL, FAS ligand; Fc, crystallizable fragment; GMP, good manufacturing practice; HDAC, histone deacetylase; hetlL-15, heterodimeric IL-15—sIL-15Ra complex; HSCT, haematopoietic stem cell transplantation; IL, interleukin; IL-2R, IL-2 receptor; IMiD, immunomodulatory drug; KIR, killer cell immunoglobulin-like receptor; mAb, monoclonal antibody; mbIL-15, membranebound IL-15; MIC, MHC class I polypeptide-related sequence; MHC, major histocompatibility complex; NAM, nicotinamide; NK, natural killer; NKG2, NK group 2; Pl, proteasome inhibitor; PD1, programmed cell death protein 1; PDL1, PD1 ligand 1; PSGL1, P-selectin glycoprotein ligand 1; RCC, renal cell carcinoma

rlL-2, recombinant IL-2; SCC, squamous cell carcinoma; TRAIL, tumour necrosis

Evidence that NK cells can

2002

mediate graft-versus-

leukaemia immunity<sup>11</sup>

factor-related apoptosis-inducing ligand, TRAILR, TRAIL receptor; T<sub>reg</sub> cell; regulatory T cell; TriKE, trispecific killer engager; scIL-15, single chain recombinant IL-15; sIL-15R $\alpha$ , soluble IL-15R $\alpha$ ; ULBP, U16-binding protein. \*Only trials studying the effect of drug treatment on NK cells as a primary or secondary end point are listed. Data presented in the tables are from ClinicalTrials.gov. †Phase II or higher.

resting NK cells14

Drugs that sensitize tumours to

TRAIL are shown to potentiate

2006

the antitumour effects of

autologous NK cells15

10. lmai, K. et al. Lancet 356, 1795–1799 (2000).

2005

1. Childs, R. W. & Carlsten, M. Nat. Rev. Drug Discov. 14, 487-498 (2015). 2. Kiessling, R. et al. Eur. J. Immunol. 5, 117–121 (1975). 3. Herberman, R. B. et al. Int. J. Cancer 16, 216–229 (1975). Koide Y. & Takasugi M. J. Natl Cancer Inst. 59, 1099–1105 (1977).
 Kärre, K. in Mechanisms of Cytotoxicity by NK Cells (eds Callewaert, D. & Herberman, R. B.) 81–91 (Academic Press, 1985). 6. Chambers W. H. et al. J. Exp. Med. 169, 1373-1389 (1989). 7. Wagtmann, N. et al. Immunity 2, 439–449 (1995). 8. Moretta, A. et al. J. Exp. Med. **182**, 875–884 (1995). 9. Braud VM et al. Nature 391, 795-799 (1998).

11. Ruggeri, L. et al. Science **295**, 2097–2100 (2002). 12. Miller, J. S. et al. Blood **105**, 3051–3057 (2005). 13. Anfossi, N. et al. Immunity 25, 331–342 (2006). 14. Bryceson, Y. T. et al. Blood 107, 159–166 (2006). 15. Lundqvist, A. et al. Cancer Res. 66, 7317–7325 (2006). 16. Berg, M. et al. Cytotherapy 11, 341-355 (2009). 17. Foley, B. et al. J. Immunol. 189, 5082-5088 (2012). 18. Carlsten, M. et al. Front. Immunol. 7, 105 (2016).

2012

Acknowledgements
This poster is adapted from the following review: Therapeutic approaches to enhance natural killer cell cytotoxicity against cancer: the force awakens.

Nat. Rev. Drug Discov. 14, 487–498 (2015). The poster content is editorially independent and the sole responsibility of Nature Publishing Group.

Edited by Alexandra Flemming; copyedited by Sophie McCabe.

© 2016 Nature Publishing Group.

http://www.nature.com/nrd/posters/nkcancertherapy

mRNA transfection of NK cells

ADCC and tumour homing<sup>18</sup>

context of HSCT<sup>1</sup>

established as a highly efficient

method to enhance the capacity of

Numerous ongoing clinical trials of

studies on the role for NK cells in the

2016

adoptive NK cell infusions and

ex vivo expanded NK cells to mediate

	Drug	Effects on NK cells	Patient populations	Comments
	es with drugs tho	nt can bolster NK ce	ell antitumour immunity*	
Cytokines	IL-2	Cytotoxicity Persistence and expansion	Melanoma, RCC, AML, neuroblastoma, breast cancer, ovarian carcinoma, fallopian tube cancer and peritoneal cancer	Some studies combine IL-2 with antitumour mAbs. rIL-2 (aldesleukin) is FDA approved
	IL-15	↑ Cytotoxicity ↑ Persistence and expansion	Melanoma, RCC, lung cancer, SCC and multiple myeloma	scIL-15 and hetIL-15 used <sup>‡</sup>
	IL-12	↑ Cytotoxicity ↑ Migration	Healthy volunteers	Lower doses improve toxicity profile
IMiDs	Lenalidomide	† Cytotoxicity † Persistence and expansion	Multiple myeloma, BCL and neuroblastoma	FDA approved
Checkpoint inhibitors	PD1-specific mAbs	↑ Cytotoxicity	Solid tumours and multiple myeloma	Tested in combination with IPH2102 (lirilumab)‡
	KIR-specific mAbs	↑ Cytotoxicity	Multiple myeloma, AML, melanoma, lung cancer and peritoneal cancer	IPH2101 and IPH2102 <sup>‡</sup>
Tumour- targeting mAbs	CD20-specific mAbs	↑ Cytotoxicity	BCL and multiple myeloma	Rituximab and veltuzumab. FDA approved
	GD2-specific mAbs	↑ Cytotoxicity	Neuroblastoma	Several different GD2-specific mAbs are being evaluated <sup>‡</sup>
	EGFR-specific mAbs	↑ Cytotoxicity	SCC	Cetuximab used in all studies <sup>‡</sup>
	ERBB2- specific mAbs	†Cytotoxicity	Breast cancer	Trastuzumab use in all studies <sup>‡</sup>
Tumour- sensitizing agents	Bortezomib	↑ Cytotoxicity	CLL, RCC, lung cancer, multiple myeloma and sarcoma	Administered before infusion of expanded NK cells to sensitize tumours to NK cell TRAIL <sup>‡</sup>
T <sub>reg</sub> cell eradication	Diphtheria toxin IL-2	↑ Cytotoxicity ↑ Persistence and expansion	AML, non-Hodgkin lymphoma and CLL	Used before NK cell infusion and in one study combined with pentostatin and rituximab <sup>‡</sup>
Clinical studi	es evaluating ad	optively infused Nk	C cells	
Activated, non- expanded NK cells	Autologous NK cells plus IL-2	↑ Cytotoxicity	Melanoma, RCC, lung cancer, nasopharyngeal cancer	Limited number studies in patien with different tumour types <sup>‡</sup>
	Autologous NK cells plus IL-15	↑ Cytotoxicity	Neuroblastoma, sarcoma, Wilms tumour and rhabdomyosarcoma	Intended to bolster NK cell tumour immunit more specifically than IL-2 does
	Allogeneic NK	↑ Cytotoxicity	AML, multiple myeloma,	Most data
	cells plus IL-2		myelodysplastic syndromes, lymphoma, ovarian carcinoma, melanoma, neuroblastoma, Ewing sarcoma, breast cancer and fallopian tube cancer	published on adoptive NK cell therapy comes from these studies <sup>‡</sup>
		↑ Cytotoxicity	myelodysplastic syndromes, lymphoma, ovarian carcinoma, melanoma, neuroblastoma, Ewing sarcoma, breast cancer and fallopian	adoptive NK cell therapy comes from these
Ex vivo expanded NK cells	Allogeneic NK cells plus	↑ Cytotoxicity  ↑ NK dose and cytotoxicity	myelodysplastic syndromes, lymphoma, ovarian carcinoma, melanoma, neuroblastoma, Ewing sarcoma, breast cancer and fallopian tube cancer	adoptive NK cell therapy comes from these studies <sup>‡</sup> Intended to bolster NK cell tumour immunit more specifically
expanded	Allogeneic NK cells plus IL-15	↑NK dose and	myelodysplastic syndromes, lymphoma, ovarian carcinoma, melanoma, neuroblastoma, Ewing sarcoma, breast cancer and fallopian tube cancer  AML and myelodysplastic syndromes  CLL, RCC, lung cancer, multiple myeloma, sarcoma, colon cancer, melanoma, neuroblastoma, prostate cancer, ALL and	adoptive NK cell therapy comes from these studies <sup>‡</sup> Intended to bolster NK cell tumour immunit more specifically than IL-2 does  Various expansion methods used, including EBV–LCL and membrane-bour cytokine/4-1BBI
expanded	Allogeneic NK cells plus IL-15  Autologous NK cells	↑NK dose and cytotoxicity  ↑NK dose and	myelodysplastic syndromes, lymphoma, ovarian carcinoma, melanoma, neuroblastoma, Ewing sarcoma, breast cancer and fallopian tube cancer  AML and myelodysplastic syndromes  CLL, RCC, lung cancer, multiple myeloma, sarcoma, colon cancer, melanoma, neuroblastoma, prostate cancer, ALL and pancreatic cancer  AML, myelodysplastic syndromes, T-cell lymphoma and	Intended to bolster NK cell tumour immunit more specifically than IL-2 does  Various expansion methods used, including EBV–LCL and membrane-bour cytokine/4-1BBI feeder cells*  Various expansion methods used, including membrane-bour cytokine/4-1BBI feeder cells. Some studuse IL-2 after